

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Cat – Domestic Short Haired

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
F0610.S9	0010076	Liver	S9	8	Mixed	1.0 mL at 20 mg/mL	400
F1501.S9	0510034	Liver	S9	1	Female	1.0 mL at 20 mg/mL	25
F1502.S9	0510084	Liver	S9	1	Female	1.0 mL at 20 mg/mL	35
F1503.S9	0510034	Liver	S9	1	Female	1.0 mL at 20 mg/mL	35
F1504.S9	0510034	Liver	S9	1	Female	1.0 mL at 20 mg/mL	80
F1001.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	35
F1002.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	45
F1003.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	30
F1004.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	35
F1000	0610343	Liver	Microsomes	3	Male	0.5 mL at 20 mg/mL	569
CP042398A	042398A	Liver	Microsomes	1	Male	0.5 mL at 20 mg/mL	9

Chicken – Broiler

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
N0610	0210299	Liver	Microsomes	22	Mixed	0.5 mL at 20 mg/mL	22
N0610	1010361	Liver	Microsomes	22	Mixed	0.5 mL at 20 mg/mL	17
N0610.S9	0210334	Liver	S9	22	Mixed	1.0 mL at 20 mg/mL	1029

Cow

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
V0610.S9	0210331	Liver	S9	2	Mixed	1.0 mL at 27 mg/mL	35
V1000.AM	PR10013	Adrenal	Mitochondria	10	Male	0.5 mL at 10 mg/mL	200
V1000.AS9	PR10014	Adrenal	S9	10	Male	1.0 mL at 10 mg/mL	400

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Dog – Beagle

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CP0710572	0710572	Intestine	S9 (No PMSF)	1	Male	1.0 mL at 4 mg/mL	52
CPPR12034	PR12034	Lung	S9 (No EDTA)	4	Mixed	1.0 mL at 5 mg/mL	95
D1000.PS	0610396	Prostate	Microsomes	5	Male	0.5 mL at 10 mg/mL	21
CP0810307	0810307	Stomach	Homogenate	3	Male	1.0 mL at 16 mg/mL	70
D1000.ZS9	PR12018	Heart	S9	2	Male	1.0 mL at 10 mg/mL	225
D1500.ZS9	PR12020	Heart	S9	2	Female	1.0 mL at 10 mg/mL	120
D1000.Z	PR12019	Heart	Microsomes	2	Male	0.5 mL at 10 mg/mL	25
D1500.Z	PR12021	Heart	Microsomes	2	Female	0.5 mL at 10 mg/mL	17
D1000.ES9	1610203	Skin	S9	3	Male	0.5 mL at 3 mg/mL	500

Gerbil – Mongolian

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
E0610	PR15038	Liver	Microsomes	60	Mixed	0.5 mL at 20 mg/mL	1
E1000.H15	PR14025	Liver	Hepatocytes	11	Male	AMY: 5.0 x 10 ⁶ per vial	50

Goat

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
C1501	013098A	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	1
C1001	061798A	Liver	Microsomes	1	Male	0.5 mL at 20 mg/mL	53
C1501.S9	013098B	Liver	S9	1	Female	1.0 mL at 20 mg/mL	7
C1502.S9	020298B	Liver	S9	1	Female	1.0 mL at 20 mg/mL	14

Guinea Pig – Hartley Albino

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
G1000.I	0510383	Intestine	Microsomes	12	Male	0.15 mL at 10 mg/mL	57
G1000.PC	0310172	Lung	Cytosol	30	Male	1.0 mL at 5 mg/mL	15
G1000.PS9	0310171	Lung	S9	30	Male	1.0 mL at 5 mg/mL	70
G1500	0910507	Liver	Microsomes	3	Female	0.5 mL at 20 mg/mL	51
G1500.S9	1510075	Liver	S9	6	Female	1.0 mL at 20 mg/mL	365
G1500.C	PR15019	Liver	Cytosol	6	Female	1.0 mL at 10 mg/mL	40

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Hamster- Golden Syrian

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
S1000.H15	1510079	Liver	Hepatocytes	6	Male	AMY: 4.0 x 10 ⁶ per vial	111
S1500.H15	PR15004	Liver	Hepatocytes	6	Female	AMY: 4.0 x 10 ⁶ per vial	128
S1500.C	PR15014	Liver	Cytosol	5	Female	1.0 mL at 10 mg/mL	80
S1500	PR15015	Liver	Microsomes	5	Female	0.5 mL at 20 mg/mL	6
S1000.ES9	PR17012	Skin	S9	8	Male	1.0 mL at 5 mg/mL	100

Horse – Quarter

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Q1501.S9	0010145	Liver	S9	1	Female	1.0 mL at 20 mg/mL	275
Q1502.S9	PR12071	Liver	S9	1	Female	1.0 mL at 20 mg/mL	49
Q1001.S9	PR12079	Liver	S9	1	Male	1.0 mL at 20 mg/mL	56
Q1501	PR09007	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	50
Q1502	PR12072	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	65
Q1001	PR12080	Liver	Microsomes	1	Male	0.5 mL at 20 mg/mL	90

Human

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
H0001.DS9	PR17027	Brain	S9	1	Male	1.0 mL at 5 mg/mL	29
H0001.DC	PR17028	Brain	Cytosol	1	Male	1.0 mL at 4 mg/mL	45
H0001.D	PR17029	Brain	Microsomes	1	Male	0.5 mL at 5 mg/mL	2
CP-16-044-Cryo	NA	Liver	Non-Parenchymal Cells	1	NA	1.0 mL	>10
H0605.H(NE)	PR15020	Liver	Homogenate (No EDTA)	10	Mixed	1.0 mL at 20 mg/mL	100
H0605.S9(NE)	1110439	Liver	S9 (No EDTA)	30	Mixed	1.0 mL at 20 mg/mL	2000
H1000.S9(NE)	PR15033	Liver	S9 (No EDTA)	4	Male	1.0 mL at 20 mg/mL	40
H1500.S9(NE)	PR15034	Liver	S9 (No EDTA)	4	Female	1.0 mL at 20 mg/mL	46
CP0710448	0710448	Liver	Microsomes (CMV-)	20	Mixed	0.5 mL at 20 mg/mL	20
CP1010169	1010169	Liver	Microsomes (Pediatric)	9	Mixed	0.5 mL at 20 mg/mL	55
CP0210296	0210296	Liver	Microsomes (High CYP3A4/5)	15	Mixed	0.5 mL at 20 mg/mL	36
CPPR09004	PR09004	Liver	Microsomes (High CYP3A4/5)	15	Mixed	0.5 mL at 20 mg/mL	25
CPPR09002	PR09002	Liver	Microsomes (High CYP3A4/5)	15	Mixed	0.5 mL at 20 mg/mL	21
CPPR12045	PR12045	Liver	Microsomes (High CYP2C19)	10	Mixed	0.5 mL at 20 mg/mL	17
CPPR13015	PR13015	Liver	Microsomes (Non-smoker)	10	Male	0.5 mL at 20 mg/mL	2
CPPR16036	PR16036	Liver	Microsomes (Non-smoker)	3	Mixed	0.5 mL at 20 mg/mL	2
CPPR16035	PR16035	Liver	Microsomes (Smoker)	3	Mixed	0.5 mL at 20 mg/mL	2

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

H0605.C(NE)	1610084	Liver	Cytosol	30	Mixed	1.0 mL at 10 mg/mL	3
H0605.RH(NE)	PR15012	Kidney	Homogenate (No EDTA)	10	Mixed	1.0 mL at 16 mg/mL	75
CP0210225	0210225	Kidney	Homogenate	4	Mixed	1.0 mL at 5 mg/mL	27
H1000.RS9	0510349	Kidney	S9	5	Male	1.0 mL at 5 mg/mL	170
H0002.R	0510087	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	2
H0004.R	0510089	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	5
H0006.R	0510091	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	1
H0008.R	0510092	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	1
H0019.R	1610057	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	9
H0020.R	1610058	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	2
H0023.R	1610059	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	6
H0024.R	1610060	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	3
H0025.R	1610061	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	13
H0026.R	1610062	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	5
H0027.R	1610063	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	5
H0028.R	1610064	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	11
H0029.R	1610065	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	9
H0030.R	1610066	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	4
H0031.R	1610067	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	3
H0032.R	1610068	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	11
CP0710571	0710571	Intestine	S9 (No PMSF)	1	Male	1.0 mL at 4 mg/mL	280
CP0910369	0910369	Intestine	Homogenate (No PMSF, Aprotinin or Leupeptin)	1	Female	1.0 mL at 4 mg/mL	350
H1000.JI	PR10076	Intestine	Microsomes (Jejunum only)	2	Male	150 µL at 10 mg/mL	103
H1000.II	PR10077	Intestine	Microsomes (Ileum only)	2	Male	150 µL at 10 mg/mL	240
H1500.JI	PR10079	Intestine	Microsomes (Jejunum only)	2	Female	150 µL at 10 mg/mL	900
H1500.II	PR10080	Intestine	Microsomes (Ileum only)	2	Female	150 µL at 10 mg/mL	365
H0093.DI	PR11017	Intestine	Microsomes (Duodenum only)	1	Female	150 µL at 10 mg/mL	1
H0093.JI	PR11018	Intestine	Microsomes (Jejunum only)	1	Female	150 µL at 10 mg/mL	65
H0093.II	PR11019	Intestine	Microsomes (Ileum only)	1	Female	150 µL at 10 mg/mL	160
H0001.LIS9	PR09019	Large Intestine	S9	1	Male	1.0 mL at 2 mg/mL	175
H0002.LIS9	0410116	Large Intestine	S9	1	Female	1.0 mL at 3 mg/mL	45
H1000.PS9(NS)	0510350	Lung	S9 (Non-Smoker)	5	Male	1.0 mL at 5 mg/mL	190
H0002.PS9	0210089	Lung	S9 (Smoker)	1	Female	1.0 mL at 5 mg/mL	65
H0005.PS9	0210091	Lung	S9 (Smoker)	1	Male	1.0 mL at 5 mg/mL	75

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

H0007.PS9	0210093	Lung	S9 (Smoker)	1	Female	1.0 mL at 5 mg/mL	100
H0008.PS9	0210095	Lung	S9 (Smoker)	1	Male	1.0 mL at 5 mg/mL	74
CPPR12028	PR12028	Lung	S9 (No EDTA, Non-Smoker)	4	Mixed	1.0 mL at 5 mg/mL	115
H0002.P	0110021	Lung	Microsomes (Smoker)	1	Female	0.5 mL at 10 mg/mL	100
H0005.P	0510090	Lung	Microsomes (Smoker)	1	Male	0.5 mL at 10 mg/mL	10
H0008.P	0210096	Lung	Microsomes (Smoker)	1	Male	0.5 mL at 10 mg/mL	10
H0001.SS9	0610086	Spleen	S9	1	Female	1.0 mL at 10 mg/mL	150
H0002.SS9	0610087	Spleen	S9	1	Male	1.0 mL at 10 mg/mL	40
H0001.ZS9	PR12022	Heart	S9	1	Male	1.0 mL at 10 mg/mL	51
H0002.ZS9	PR12024	Heart	S9	1	Female	1.0 mL at 10 mg/mL	61
H1000.U	PR16025	Muscle	Microsomes	3	Mixed	0.5 mL at 10 mg/mL	26

Minipig – Gottingen

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Z6500.H15	PR15006	Liver	Hepatocytes	1	Female	AMY: 4.0 x 10 ⁶ per vial	9
Z6511.S9	PR15016	Liver	S9	1	Female	1.0 mL at 20 mg/mL	65
Z6511.C	PR15017	Liver	Cytosol	1	Female	1.0 mL at 10 mg/mL	55
Z6511	PR15018	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	62
Z6500.YS9	PR14037	Eye	S9	10	Female	1.0 mL at 5 mg/mL	41

Minipig – Hanford

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Z5001.S9	0710107	Liver	S9	1	Male	1.0 mL at 20 mg/mL	7

Minipig – Sinclair

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Z2000.IS9	0310177	Intestine	S9	3	Male	1.0 mL at 4 mg/mL	230
Z2000.C	0310182	Liver	Cytosol	3	Male	1.0 mL at 10 mg/mL	160

Minipig – Yucatan

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Z3000.IS9	0310178	Intestine	S9	3	Male	1.0 mL at 4 mg/mL	166
Z3000.C	0310185	Liver	Cytosol	3	Male	1.0 mL at 10 mg/mL	190

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Monkey – Rhesus

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
P1000.IC	0410224	Intestine	Cytosol	7	Male	1.0 mL at 3 mg/mL	35
P1000.I	0410225	Intestine	Microsomes	7	Male	0.15 mL at 10 mg/mL	3
P1000.I	1610411	Intestine	Microsomes	7	Male	0.15 mL at 10 mg/mL	62
P1000.IS9	0410138	Intestine	S9	7	Male	1.0 mL at 4 mg/mL	210
P1000.PC	0310160	Lung	Cytosol	4	Male	1.0 mL at 5 mg/mL	20
P1000.P	0510272	Lung	Microsomes	4	Male	0.5 mL at 5 mg/mL	15
P1000.RS9	0110170	Kidney	S9	2	Male	1.0 mL at 5 mg/mL	18

Monkey – Cynomolgus

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR12033	PR12033	Lung	S9 (No EDTA)	4	Mixed	1.0 mL at 5 mg/mL	100
P2012.RS9	PR10060	Kidney	S9	1	Male	1.0 mL at 5 mg/mL	40
P2013.RS9	PR10061	Kidney	S9	1	Male	1.0 mL at 5 mg/mL	40
P2034.RS9	PR10062	Kidney	S9	1	Male	1.0 mL at 5 mg/mL	30
P2000.RH	PR14026	Kidney	Homogenate	10	Male	1.0 mL at 5 mg/mL	2
CP0610033	0610033	Spleen	S9	2	Male	1.0 mL at 10 mg/mL	17
CP0610068	0610068	Brain	Mitochondria	1	Male	0.5 mL at 10 mg/mL	20
P2500.S9(NE)	PR15032	Liver	S9 (No EDTA)	3	Female	1.0 mL at 20 mg/mL	23
P2000.AS9	PR16034	Adrenal	S9	10	Male	1.0 mL at 5 mg/mL	30

Mouse – BALB/c

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M3500.H15	1210157	Liver	Hepatocytes	Pooled	Female	AMY: 2.0 x 10 ⁶ per vial	2
M3500.H15	1610338	Liver	Hepatocytes	Pooled	Female	AMY: 2.0 x 10 ⁶ per vial	8
M3500.PS9	0610103	Lung	S9	26	Female	1.0 mL at 5 mg/mL	15
M3000.IS9	PR13010	Intestine	S9	60	Male	1.0 mL at 4 mg/mL	18

Mouse – B6C3F1

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M2000.H15	1510089	Liver	Hepatocytes	24	Male	AMY: 2.0 x 10 ⁶ per vial	55
M2500.H15	1510123	Liver	Hepatocytes	24	Female	AMY: 2.0 x 10 ⁶ per vial	13
M2500.C	PR14038	Liver	Cytosol	30	Female	1.0 mL at 10 mg/mL	48
M2500	PR14039	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	13

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Mouse – B6SJLF1

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CP0710313	0710313	Liver	Microsomes	146	Male	0.5 mL at 20 mg/mL	177

Mouse – CD1

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M1114	0710616	Liver	Microsomes (Phenobarbital Treated)	50	Male	0.5 mL at 20 mg/mL	14
M1114.S9	0710618	Liver	S9 (Phenobarbital Treated)	50	Male	1.0 mL at 20 mg/mL	113
M0610	PR10002	Liver	Microsomes	1609	Mixed	0.5 mL at 20 mg/mL	2
M1000.PH	0110114	Lung	Homogenate	196	Male	1.0 mL at 5 mg/mL	14
M1500.PS9	PR10084	Lung	S9	101	Female	1.0 mL at 5 mg/mL	44
CPPR12026	PR12026	Lung	S9 (No EDTA)	30	Mixed	1.0 mL at 5 mg/mL	55
M1500.RS9	PR10085	Kidney	S9	101	Female	1.0 mL at 5 mg/mL	400
CPPR11006	PR11006	Intestine	S9 (No EDTA - scraped)	41	Male	1.0 mL at 4 mg/mL	55
M1500.IS9	PR10083	Intestine	S9	101	Female	1.0 mL at 4 mg/mL	14

Mouse – C57BL/6

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M5000.S9(NE)	PR15029	Liver	S9 (No EDTA)	5	Male	1.0 mL at 20 mg/mL	3
M5000.LT	PR16037	Liver	Tritosomes	47	Male	0.25 mL at 3.3 mg/mL	2
M5000.IS9	PR13011	Intestine	S9	57	Male	1.0 mL at 4 mg/mL	20

Mouse – C57BL/6J

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M6500.H15	PR14031	Liver	Hepatocytes	8	Female	AMY: 2.0 x 10 ⁶ per vial	2
M6500.PS9	PR14034	Lung	S9	20	Female	1.0 mL at 5 mg/mL	24

Mouse – db/db Obese

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09048	PR09048	Liver	Microsomes	20	Male	0.5 mL at 20 mg/mL	28
CPPR09050	PR09050	Liver	Microsomes	20	Female	0.5 mL at 20 mg/mL	2

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Mouse – Fox Chase SCID

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CP0510352	0510352	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	6
CP0810346	0810346	Liver	Microsomes	60	Female	0.5 mL at 20 mg/mL	5
CPPR12064	PR12064	Liver	Microsomes	12	Mixed	0.5 mL at 20 mg/mL	4
CPPR12066	PR12066	Intestine	Microsomes	50	Female	0.15 mL at 10 mg/mL	3

Mouse – KK.Cg A^v/J

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09040	PR09040	Liver	Microsomes	20	Male	0.5 mL at 20 mg/mL	16
CPPR09042	PR09042	Liver	Microsomes	20	Female	0.5 mL at 20 mg/mL	25

Mouse – KK/HIJ

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09041	PR09041	Liver	Microsomes	20	Male	0.5 mL at 20 mg/mL	26
CPPR09043	PR09043	Liver	Microsomes	20	Female	0.5 mL at 20 mg/mL	16

Mouse – ob/ob Lean

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09045	PR09045	Liver	Microsomes	20	Male	0.5 mL at 20 mg/mL	3
CPPR09057	PR09057	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	2

Mouse – ob/ob Obese

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09044	PR09044	Liver	Microsomes	20	Male	0.5 mL at 20 mg/mL	26
CPPR09046	PR09046	Liver	Microsomes	20	Female	0.5 mL at 20 mg/mL	21

Mouse – Swiss Webster

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CP0510265	0510265	Brain	S9	40	Male	1.0 mL at 10 mg/mL	44
M4000	0810446	Liver	Microsomes	150	Male	0.5 mL at 20 mg/mL	150

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Pig – Standard

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Z1503	060998A	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	17
Z1501	011498A	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	28
Z0610	0210297	Liver	Microsomes	2	Mixed	0.5 mL at 20 mg/mL	56
Z1001.S9	011498B	Liver	S9	1	Male	1.0 mL at 20 mg/mL	170
Z0610.S9	0210330	Liver	S9	2	Mixed	1.0 mL at 20 mg/mL	430

Rabbit – Dutch Belted

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
L2500	0410205	Liver	Microsomes	3	Female	0.5 mL at 20 mg/mL	34
L2500.S9	0710631	Liver	S9	3	Female	1.0 mL at 20 mg/mL	244

Rabbit – New Zealand

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
L1500.I	PR10041	Intestine	Microsomes	2	Female	0.15 mL at 10 mg/mL	9
L1000.ZS9	PR12014	Heart	S9	4	Male	1.0 mL at 10 mg/mL	1
L1500.ZS9	PR12016	Heart	S9	4	Female	1.0 mL at 10 mg/mL	3
L1000.Z	PR12015	Heart	Microsomes	4	Male	0.5 mL at 10 mg/mL	3
L1500.Z	PR12017	Heart	Microsomes	4	Female	0.5 mL at 10 mg/mL	6
L1000.YS9	PR12052	Eye	S9	75	Male	0.5 mL at 5 mg/mL	530

Rat – Fischer

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
R2000.RS9	0710638	Kidney	S9	9	Male	1.0 mL at 5 mg/mL	75
R2000.IS9	0710637	Intestine	S9	10	Male	1.0 mL at 5 mg/mL	24
R2000.P	0810421	Lung	Microsomes	30	Male	0.5 mL at 10 mg/mL	8

Rat – GK

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09036	PR09036	Liver	Microsomes	3	Male	0.5 mL at 20 mg/mL	18

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Rat – Lewis

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
R5000.S9	0410316	Liver	S9	3	Male	1.0 mL at 20 mg/mL	4
R5000	0910114	Liver	Microsomes	3	Male	0.5 mL at 20 mg/mL	1
R5500	PR10086	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	207

Rat – Long Evans

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CP0210409	0210409	Liver	Microsomes	10	Male	0.5 mL at 20 mg/mL	130

Rat – Sprague-Dawley

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR11007	PR11007	Liver	S9 (No EDTA)	10	Male	1.0 mL at 20 mg/mL	16
R1500.S9(NE)	PR15031	Liver	S9 (No EDTA)	5	Female	1.0 mL at 20 mg/mL	145
R1000.M	1010131	Liver	Mitochondria	5	Male	0.5 mL at 20 mg/mL	3
R1000.M	1110056	Liver	Mitochondria	15	Male	0.5 mL at 20 mg/mL	97
R0610	PR10001	Liver	Microsomes	515	Mixed	1.0 mL at 20 mg/mL	2
R0605.C(NE)	PR16028	Liver	Cytosol	15	Mixed	1.0 mL at 10 mg/mL	3
R1000.L	1710154	Liver	Lysosomes	4	Male	0.25 mL at 1.7 mg/mL	51
R1000.DH(NE)	1810159	Brain	Homogenate	12	Male	1.0 mL at 5 mg/mL	30
CP0510266	0510266	Brain	S9	30	Male	1.0 mL at 10 mg/mL	100
CP0610001	0610001	Testicle	Microsomes	250	Male	0.5 mL at 20 mg/mL	3
R1000.THO	PR12035	Testicle	Homogenate	75	Male	0.5 mL at 10 mg/mL	160
R1000.TS9	PR12036	Testicle	S9	75	Male	0.5 mL at 10 mg/mL	1060
R1000.PH	0710553	Lung	Homogenate	25	Male	1.0 mL at 20 mg/mL	5
R1000.PH	0110118	Lung	Homogenate	24	Male	1.0 mL at 5 mg/mL	11
CPPR12027	PR12027	Lung	S9 (No EDTA)	10	Mixed	1.0 mL at 5 mg/mL	140
CPPR11093	PR11093	Intestine	S9 (No PMSF)	40	Male	1.0 mL at 4 mg/mL	59
R0610.ZM	PR10055	Heart	Mitochondria	250	Mixed	0.5 mL at 20 mg/mL	414
R0610.ZS9	PR10056	Heart	S9	250	Mixed	1.0 mL at 20 mg/mL	390
R1500.ZS9	PR12008	Heart	S9	60	Female	1.0 mL at 10 mg/mL	2
R1500.Z	PR12009	Heart	Microsomes	60	Female	0.5 mL at 10 mg/mL	3
R1000.YS9	PR12053	Eye	S9	200	Male	0.5 mL at 5 mg/mL	620
CP0610074	0610074	Adrenal	Microsomes	750	Female	0.5 mL at 10 mg/mL	2

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Rat – Sprague-Dawley - Treated

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
R1058	0494	Liver	Microsomes (PFDA-treated)	15	Male	0.5 mL at 10 mg/mL	235
R1063.S9	0210128	Liver	S9 (Clofibrac acid-treated)	25	Male	1.0 mL at 20 mg/mL	200
R1078.S9	0110142	Liver	S9 (Phenobarbital-treated)	24	Male	1.0 mL at 20 mg/mL	140
R1084	0494	Liver	Microsomes (3MC-treated)	16	Male	0.5 mL at 10 mg/mL	2
R1084.S9	0210131	Liver	S9 (3MC-treated)	25	Male	1.0 mL at 20 mg/mL	800
R1088.S9	0210134	Liver	S9 (Isoniazid-treated)	25	Male	1.0 mL at 20 mg/mL	60
R1073.DH	PR17014	Brain	Homogenate (Saline-treated)	12	Male	1.0 mL at 5 mg/mL	4
R1078.DH	PR17013	Brain	Homogenate (Phenobarbital-treated)	12	Male	1.0 mL at 5 mg/mL	4
R1073.DS9	PR17016	Brain	S9 (Saline-treated)	12	Male	1.0 mL at 5 mg/mL	4
R1078.DS9	PR17015	Brain	S9 (Phenobarbital-treated)	12	Male	1.0 mL at 5 mg/mL	4
R1073.D	PR17018	Brain	Microsomes (Saline-treated)	12	Male	0.5 mL at 5 mg/mL	2
R1078.D	PR17017	Brain	Microsomes (Phenobarbital-treated)	12	Male	0.5 mL at 5 mg/mL	2

Rat – Wistar

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
R3500.H15	PR14042	Liver	Hepatocytes	7	Female	AMY: 7.0 x 10 ⁶ per vial	28
R3500.H15	PR15021	Liver	Hepatocytes	10	Female	AMY: 7.0 x 10 ⁶ per vial	53
CP-14-055	PR14040	Liver	Hepatocytes (Harlan)	8	Male	AMY: 7.0 x 10 ⁶ per vial	47

Rat – Wistar Han

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR13033	PR13033	Intestine	Microsomes (No PMSF)	75	Male	0.15 mL at 10 mg/mL	108
R6000.IC(NP)	PR16026	Intestine	Cytosol (No PMSF)	10	Male	1.0 mL at 4 mg/mL	14
CPPR17023	PR17023	Kidney	Membrane	6	Male	0.7 mL at 22 mg/mL	1
R6500	1410203	Liver	Microsomes	225	Female	0.5 mL at 20 mg/mL	950
CPPR16054	PR16054	Liver	Microsomes	40	Mixed	0.5 mL at 20 mg/mL	370
CPPR16053	PR16053	Liver	S9	40	Mixed	1.0 mL at 20 mg/mL	52

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 29 JUNE 2018

Rat – ZDF Lean

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09053	PR09053	Liver	Microsomes	6	Male	0.5 mL at 20 mg/mL	8
CPPR09055	PR09055	Liver	Microsomes	7	Female	0.5 mL at 20 mg/mL	4

Rat – ZDF Obese

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR09032	PR09032	Liver	Microsomes	3	Male	0.5 mL at 20 mg/mL	4
CPPR09054	PR09054	Liver	Microsomes	5	Female	0.5 mL at 20 mg/mL	5

Sheep

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
O1000.S9	0410144	Liver	S9	2	Male	1.0 mL at 20 mg/mL	147
O1000.P	0210188	Lung	Microsomes	2	Male	0.5 mL at 10 mg/mL	22
O1000.PS9	0210187	Lung	S9	2	Male	1.0 mL at 5 mg/mL	195

These products are surplus stock of previously prepared custom products. These products are not included in Sekisui XenoTech's standard product line. As such, their inventory stock is not renewed. Please contact Customer Service to confirm availability. Future orders of the above listed products may incur custom product processing fees if new product preparation is required. Please allow up to 5 days for dispensing and packaging.