

## Surplus Custom Products

**AVAILABLE AS OF 17 MARCH 2021**

---

### 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	0510034	Liver	S9	1	Female	1.0 mL at 20 mg/mL	30
F1503.S9	0510034	Liver	S9	1	Female	1.0 mL at 20 mg/mL	25
F1504.S9	0510034	Liver	S9	1	Female	1.0 mL at 20 mg/mL	65
F1001.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	25
F1002.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	40
F1003.S9	0510034	Liver	S9	1	Male	1.0 mL at 20 mg/mL	25
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	350

### Chicken – Broiler

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
N0610	1910153	Liver	Microsomes	22	Mixed	0.5 mL at 20 mg/mL	130
N0610.S9	0210334	Liver	S9	22	Mixed	1.0 mL at 20 mg/mL	440

### 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	14
V1000.AM	PR10013	Adrenal	Mitochondria	10	Male	0.5 mL at 10 mg/mL	150
V1000.AS9	PR10014	Adrenal	S9	10	Male	1.0 mL at 10 mg/mL	400

### Dog – Beagle

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR12034	PR12034	Lung	S9 (No EDTA)	4	Mixed	1.0 mL at 5 mg/mL	92
D1000.ZS9	PR12018	Heart	S9	2	Male	1.0 mL at 10 mg/mL	200
D1500.ZS9	PR12020	Heart	S9	2	Female	1.0 mL at 10 mg/mL	110
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	1.0 mL at 3 mg/mL	188

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 17 MARCH 2021**

### Gerbil – Mongolian

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
E1000.H15	PR14025	Liver	Hepatocytes	11	Male	AMY: 5.0 x 10 <sup>6</sup> 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	4
C1502.S9	020298B	Liver	S9	1	Female	1.0 mL at 20 mg/mL	4

### 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	5
G1000.PS9	0310171	Lung	S9	30	Male	1.0 mL at 5 mg/mL	50
G1500	0910507	Liver	Microsomes	3	Female	0.5 mL at 20 mg/mL	36
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	70

### 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 <sup>6</sup> per vial	107
S1500.H15	PR15004	Liver	Hepatocytes	6	Female	AMY: 4.0 x 10 <sup>6</sup> per vial	125
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	155

### 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	273
Q1502.S9	PR12071	Liver	S9	1	Female	1.0 mL at 20 mg/mL	47
Q1001.S9	PR12079	Liver	S9	1	Male	1.0 mL at 20 mg/mL	55
Q1502	PR12072	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	7
Q1001	PR12080	Liver	Microsomes	1	Male	0.5 mL at 20 mg/mL	90

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 17 MARCH 2021**

### 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	9
H0001.DC	PR17028	Brain	Cytosol	1	Male	1.0 mL at 4 mg/mL	23
H0605.H(NE)	PR15020	Liver	Homogenate (No EDTA)	10	Mixed	1.0 mL at 20 mg/mL	27
CPPR20016	PR20016	Liver	Homogenate- custom buffer	6	Mixed	1.0 mL	5
CPPR20019	PR20019	Liver	Crude Homogenate	3	Male	1.0 mL at 20 mg/mL	17
H0605.S9(NE)	1110439	Liver	S9 (No EDTA)	30	Mixed	1.0 mL at 20 mg/mL	1968
H1500.S9(NE)	PR15034	Liver	S9 (No EDTA)	4	Female	1.0 mL at 20 mg/mL	46
CP1010169	1010169	Liver	Microsomes (Pediatric)	9	Mixed	0.5 mL at 20 mg/mL	55
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
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	58
CP0210225	0210225	Kidney	Homogenate	4	Mixed	1.0 mL at 5 mg/mL	19
H1000.RS9	0510349	Kidney	S9	5	Male	1.0 mL at 5 mg/mL	170
H0004.R	0510089	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	2
H0019.R	1610057	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	7
H0023.R	1610059	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	4
H0024.R	1610060	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	1
H0025.R	1610061	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	11
H0026.R	1610062	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	3
H0027.R	1610063	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	3
H0028.R	1610064	Kidney	Microsomes	1	Male	0.5 mL at 10 mg/mL	8
H0029.R	1610065	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	7
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	1
H0032.R	1610068	Kidney	Microsomes	1	Female	0.5 mL at 10 mg/mL	8
H0610.IH	PR14023	Intestine	Homogenate	12	Mixed	1.0 mL at 4 mg/mL	33
CP0710571	0710571	Intestine	S9 (No PMSF)	1	Male	1.0 mL at 4 mg/mL	276
H1000.JI	PR10076	Intestine	Microsomes (Jejunum only)	2	Male	150 µL at 10 mg/mL	102
H1000.II	PR10077	Intestine	Microsomes (Ileum only)	2	Male	150 µL at 10 mg/mL	237
H1500.JI	PR10079	Intestine	Microsomes (Jejunum only)	2	Female	150 µL at 10 mg/mL	899
H1500.II	PR10080	Intestine	Microsomes (Ileum only)	2	Female	150 µL at 10 mg/mL	362

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 17 MARCH 2021**

H0093.DI	PR11017	Intestine	Microsomes (Duodenum only)	1	Female	150 µL at 10 mg/mL	8
H0093.JI	PR11018	Intestine	Microsomes (Jejunum only)	1	Female	150 µL at 10 mg/mL	64
H0093.II	PR11019	Intestine	Microsomes (Ileum only)	1	Female	150 µL at 10 mg/mL	120
H0001.LIS9	PR09019	Large Intestine	S9	1	Male	1.0 mL at 2 mg/mL	124
H0002.LIS9	0410116	Large Intestine	S9	1	Female	1.0 mL at 3 mg/mL	43
H1000.PS9(NS)	0510350	Lung	S9 (Non-Smoker)	5	Male	1.0 mL at 5 mg/mL	180
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
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	112
H0610.QS9	2010063	Pancreas	S9	2	Mixed	1.0 mL at 4 mg/mL	48
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	38
H0001.ZS9	PR12022	Heart	S9	1	Male	1.0 mL at 10 mg/mL	42
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	19

### Minipig – Gottingen

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
Z6004.IS9(NP)	2010222	Intestine	S9- No PMSF	1	Male	1.0 mL at 4 mg/mL	42
Z6000.RS9	2010134	Kidney	S9	3	Male	1.0 mL at 5 mg/mL	48
Z6500.H15	PR15006	Liver	Hepatocytes	1	Female	AMY: 4.0 x 10 <sup>6</sup> per vial	5
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 – 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	17
Z2000.C	0310182	Liver	Cytosol	3	Male	1.0 mL at 10 mg/mL	130

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 17 MARCH 2021**

### 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	140
Z3000.C	0310185	Liver	Cytosol	3	Male	1.0 mL at 10 mg/mL	190

### 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	50
P1000.IS9	0410138	Intestine	S9	7	Male	1.0 mL at 4 mg/mL	280
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
P2610.S9(NE)	PR19032	Liver	S9 (No EDTA)	2	Mixed	1.0 mL at 20 mg/mL	12
P2006A.L	1810033	Liver	Lysosomes	1	Male	0.25 mL at 6 mg/mL	10
P2000.M	2010229	Liver	Mitochondria	3	Male	0.5 mL at 20 mg/mL	38
P2610.M	2010227	Liver	Mitochondria	5	Mixed	1.0 mL at 20 mg/mL	3
P2610.S9	2010226	Liver	S9	5	Mixed	1.0 mL at 17 mg/mL	58
CPPR12033	PR12033	Lung	S9 (No EDTA)	4	Mixed	1.0 mL at 5 mg/mL	94
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	30
P2000.AS9	PR16034	Adrenal	S9	10	Male	1.0 mL at 5 mg/mL	30
P0610.QS9	2010100	Pancreas	S9	2	Mixed	1.0 mL at 4 mg/mL	75

### 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 <sup>6</sup> per vial	2
M3500.H15	1610338	Liver	Hepatocytes	Pooled	Female	AMY: 2.0 x 10 <sup>6</sup> per vial	7
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

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 17 MARCH 2021**

### 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 <sup>6</sup> per vial	53
M2500.H15	1510123	Liver	Hepatocytes	24	Female	AMY: 2.0 x 10 <sup>6</sup> per vial	11
M2500.C	PR14038	Liver	Cytosol	30	Female	1.0 mL at 10 mg/mL	47
M2500	PR14039	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	12

### 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	60
M1114.S9	0710618	Liver	S9 (Phenobarbital Treated)	50	Male	1.0 mL at 20 mg/mL	150
M0610.S9	2010224	Liver	S9	15	Mixed	1.0 mL at 17 mg/mL	55
M1000.M	2010230	Liver	Mitochondria	30	Male	0.5 mL at 20 mg/mL	125
M1500.PS9	PR10084	Lung	S9	101	Female	1.0 mL at 5 mg/mL	14
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	397
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	10
M0610.QS9	2010061	Pancreas	S9	50	Mixed	1.0 mL at 4 mg/mL	162

### Mouse – C57BL/6

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M5000.IS9	PR13011	Intestine	S9	57	Male	1.0 mL at 4 mg/mL	20
M5610.M	2010225	Liver	Mitochondria	15	Mixed	1.0 mL at 20 mg/mL	17
M5000.L	1810068	Liver	Lysosomes	38	Male	0.25 mL at 1.8 mg/mL	49

### Mouse – C57BL/6J

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
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 17 MARCH 2021**

---

### Mouse – Fox Chase SCID

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR12066	PR12066	Intestine	Microsomes	50	Female	0.15 mL at 10 mg/mL	3

### Mouse – KK.Cg A<sup>y</sup>/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	15
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	25
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	2
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	24
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	31
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 17 MARCH 2021

### 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	47
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	425

### 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	6
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	430

### 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	70
R2000.P	0810421	Lung	Microsomes	30	Male	0.5 mL at 10 mg/mL	8

### Rat – Lewis

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
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	194

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 17 MARCH 2021**

### 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	65
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	65
CPPR20020	PR20020	Liver	Crude Homogenate	3	Male	1.0 mL at 20 mg/mL	15
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	50
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	26
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	150
R1000.TS9	PR12036	Testicle	S9	75	Male	0.5 mL at 10 mg/mL	1055
CPPR12027	PR12027	Lung	S9 (No EDTA)	10	Mixed	1.0 mL at 5 mg/mL	137
R1000.PH	2010231	Lung	Homogenate	25	Male	1.0 mL at 12 mg/mL	86
R1000.PH	2010232	Lung	Homogenate	25	Male	1.0 mL at 12 mg/mL	24
CPPR11093	PR11093	Intestine	S9 (No PMSF)	40	Male	1.0 mL at 4 mg/mL	55
R0610.QS9	2010062	Pancreas	S9	4	Mixed	1.0 mL at 4 mg/mL	18
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	380
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	597
CPPR19020	PR19020	Kidney	S9 (No EDTA)	8	Male	1.0 mL at 5 mg/mL	100
R1000.RH	2010233	Kidney	Homogenate	25	Male	1.0 mL at 20 mg/mL	180

### 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	140
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

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 17 MARCH 2021**

R1078.DH	PR17013	Brain	Homogenate (Phenobarbital-treated)	12	Male	1.0 mL at 5 mg/mL	4
----------	---------	-------	---------------------------------------	----	------	-------------------	---

**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 <sup>6</sup> per vial	23
R3500.H15	PR15021	Liver	Hepatocytes	10	Female	AMY: 7.0 x 10 <sup>6</sup> per vial	21
CP-14-055	PR14040	Liver	Hepatocytes (Harlan)	8	Male	AMY: 7.0 x 10 <sup>6</sup> per vial	47

**Rat – Wistar Han**

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
CPPR20014	PR20014	Intestine	S9 (No PMSF)	25	Male	1.0 mL at 4 mg/mL	41
CPPR13033	PR13033	Intestine	Microsomes (No PMSF)	75	Male	0.15 mL at 10 mg/mL	97
R6000.IC(NP)	PR16026	Intestine	Cytosol (No PMSF)	10	Male	1.0 mL at 4 mg/mL	14
R6500	1410203	Liver	Microsomes	225	Female	0.5 mL at 20 mg/mL	940

**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	7
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
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	105
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.