

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

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	42
F1000	0610343	Liver	Microsomes	3	Male	0.5 mL at 20 mg/mL	315

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	150
N0610.S9	0210334	Liver	S9	22	Mixed	1.0 mL at 20 mg/mL	565

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.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	26
D1000.YS9	2110297	Eye	S9	3	Male	0.5 mL at 5 mg/mL	35
D1000.ZS9	PR12018	Heart	S9	2	Male	1.0 mL at 10 mg/mL	12
D1500.ZS9	PR12020	Heart	S9	2	Female	1.0 mL at 10 mg/mL	20
D1000.Z	PR12019	Heart	Microsomes	2	Male	0.5 mL at 10 mg/mL	30
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	190

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 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	4
C1502.S9	020298B	Liver	S9	1	Female	1.0 mL at 20 mg/mL	4

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

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	20
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	39
G1500.S9	1510075	Liver	S9	6	Female	1.0 mL at 20 mg/mL	20
G1500.C	PR15019	Liver	Cytosol	6	Female	1.0 mL at 10 mg/mL	40

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	37
S1500.H15	PR15004	Liver	Hepatocytes	6	Female	AMY: 4.0 x 10 ⁶ per vial	45
S1500	PR21017	Liver	Microsomes	6	Female	0.5 mL at 20 mg/mL	27
S1000.ES9	PR17012	Skin	S9	8	Male	1.0 mL at 5 mg/mL	40

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	20
Q1001.S9	PR12079	Liver	S9	1	Male	1.0 mL at 20 mg/mL	1
Q1001	PR12080	Liver	Microsomes	1	Male	0.5 mL at 20 mg/mL	37

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

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	2
H0001.DC	PR17028	Brain	Cytosol	1	Male	1.0 mL at 4 mg/mL	17
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	217
CP1010169	1010169	Liver	Microsomes (Pediatric)	9	Mixed	0.5 mL at 20 mg/mL	57
CPPR12045	PR12045	Liver	Microsomes (High CYP2C19)	10	Mixed	0.5 mL at 20 mg/mL	17
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	18
CP0210225	0210225	Kidney	Homogenate	4	Mixed	1.0 mL at 5 mg/mL	9
H1000.RS9	0510349	Kidney	S9	5	Male	1.0 mL at 5 mg/mL	165
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	6
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	9
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	2
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	28
H1000.PS9(NS)	0510350	Lung	S9 (Non-Smoker)	5	Male	1.0 mL at 5 mg/mL	80
CPPR12028	PR12028	Lung	S9 (No EDTA, Non-Smoker)	4	Mixed	1.0 mL at 5 mg/mL	106
H0610.QS9	2010063	Pancreas	S9	2	Mixed	1.0 mL at 4 mg/mL	52
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	41
H0002.ZS9	PR12024	Heart	S9	1	Female	1.0 mL at 10 mg/mL	12
H1000.U	PR16025	Muscle	Microsomes	3	Mixed	0.5 mL at 10 mg/mL	19

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

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	27
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 ⁶ 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	20
Z6511	PR15018	Liver	Microsomes	1	Female	0.5 mL at 20 mg/mL	60

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

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.IS9	0410138	Intestine	S9	7	Male	1.0 mL at 4 mg/mL	30
P1000.RS9	0110170	Kidney	S9	2	Male	1.0 mL at 5 mg/mL	18
P1000	1110332	Liver	Microsomes	6	Male	0.5 mL at 20 mg/mL	480
P1000.C	1110331	Liver	Cytosol	6	Male	1.0 mL at 10 mg/mL	175
P1000.C	1810192	Liver	Cytosol	3	Male	1.0 mL at 10 mg/mL	175
P1000.S9	1110330	Liver	S9	6	Male	1.0 mL at 20 mg/mL	36
P1500	1210053	Liver	Microsomes	6	Female	0.5 mL at 20 mg/mL	403
P1500.C	0710096	Liver	Cytosol	7	Female	1.0 mL at 10 mg/mL	85
P1500.C	1210052	Liver	Cytosol	6	Female	1.0 mL at 10 mg/mL	154
P1500.S9	1110079	Liver	S9	5	Female	1.0 mL at 20 mg/mL	56

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

Monkey – Cynomolgus

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
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	23
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
CP0610033	0610033	Spleen	S9	2	Male	1.0 mL at 10 mg/mL	25
P2000.AS9	PR16034	Adrenal	S9	10	Male	1.0 mL at 5 mg/mL	28
P0610.QS9	2010100	Pancreas	S9	2	Mixed	1.0 mL at 4 mg/mL	79

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	5
M3500.H15	1610338	Liver	Hepatocytes	Pooled	Female	AMY: 2.0 x 10 ⁶ per vial	3

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	53
M2500.H15	1510123	Liver	Hepatocytes	24	Female	AMY: 2.0 x 10 ⁶ per vial	11
M2500	PR14039	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	11

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	55
M1114.S9	0710618	Liver	S9 (Phenobarbital Treated)	50	Male	1.0 mL at 20 mg/mL	70
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	10
M1500.RS9	PR10085	Kidney	S9	101	Female	1.0 mL at 5 mg/mL	30
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	166

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

Mouse – C57BL/6

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
M5610.M	2010225	Liver	Mitochondria	15	Mixed	1.0 mL at 20 mg/mL	7
M5000.L	1810068	Liver	Lysosomes	38	Male	0.25 mL at 1.8 mg/mL	300

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

Pig – Standard

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
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	3
Z0610.S9	0210330	Liver	S9	2	Mixed	1.0 mL at 20 mg/mL	55

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	3
L1000.YS9	PR12052	Eye	S9	75	Male	0.5 mL at 5 mg/mL	200

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	42

Rat – Lewis

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
R5500	PR10086	Liver	Microsomes	30	Female	0.5 mL at 20 mg/mL	20

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

Rat – Sprague-Dawley

Product ID	Lot	Organ	Subcellular Fraction	Pool size	Gender	Vial Size	Vials Available
R1500.S9(NE)	PR15031	Liver	S9 (No EDTA)	5	Female	1.0 mL at 20 mg/mL	60
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	20
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	30
R1000.DH(NE)	1810159	Brain	Homogenate	12	Male	1.0 mL at 5 mg/mL	81
CP0510266	0510266	Brain	S9	30	Male	1.0 mL at 10 mg/mL	17
CP0610001	0610001	Testicle	Microsomes	250	Male	0.5 mL at 20 mg/mL	3
R1000.TS9	PR12036	Testicle	S9	75	Male	0.5 mL at 10 mg/mL	302
CPPR12027	PR12027	Lung	S9 (No EDTA)	10	Mixed	1.0 mL at 5 mg/mL	16
R1000.PH	2010231	Lung	Homogenate	25	Male	1.0 mL at 12 mg/mL	81
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	54
R0610.QS9	2010062	Pancreas	S9	4	Mixed	1.0 mL at 4 mg/mL	18
R0610.ZS9	PR10056	Heart	S9	250	Mixed	1.0 mL at 20 mg/mL	30
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	98
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
R1063.S9	0210128	Liver	S9 (Clofibrac acid-treated)	25	Male	1.0 mL at 20 mg/mL	200
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	900
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

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	19
R3500.H15	PR15021	Liver	Hepatocytes	10	Female	AMY: 7.0 x 10 ⁶ per vial	21

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.

Surplus Custom Products

AVAILABLE AS OF 17 FEBRUARY 2023

Rat – Wistar Han

<u>Product ID</u>	<u>Lot</u>	<u>Organ</u>	<u>Subcellular Fraction</u>	<u>Pool size</u>	<u>Gender</u>	<u>Vial Size</u>	<u>Vials Available</u>
CPPR20014	PR20014	Intestine	S9 (No PMSF)	25	Male	1.0 mL at 4 mg/mL	28
CPPR13033	PR13033	Intestine	Microsomes (No PMSF)	75	Male	0.15 mL at 10 mg/mL	22
R6500	1410203	Liver	Microsomes	225	Female	0.5 mL at 20 mg/mL	1

Sheep

<u>Product ID</u>	<u>Lot</u>	<u>Organ</u>	<u>Subcellular Fraction</u>	<u>Pool size</u>	<u>Gender</u>	<u>Vial Size</u>	<u>Vials Available</u>
O1000.S9	0410144	Liver	S9	2	Male	1.0 mL at 20 mg/mL	12

These products are surplus stock of previously prepared custom products. These products are not included in 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 7 days for dispensing and packaging.