

LEXSY Products – Overview

LEXSY Expression Kits

Product	Cat. No.	Amount	Price EUR
LEXSYcon2 Expression Kit for constitutive protein expression, contains vector pLEXSY-sat2 or pLEXSY-ble2 or pLEXSY-hyg2 or pLEXSY-neo2	EGE-1300	1 Kit	960,-
LEXSInduce3 Expression Kit for inducible expression from integrative constructs, contains vector pLEXSY_I-blecherry3 or pLEXSY_I-ble3 or pLEXSY_I-neo3	EGE-1410	1 Kit	2250,-
LEXSInduce4 Expression Kit for inducible expression from episomal constructs, contains vector pLEXSY_IE-blecherry4	EGE-1420	1 Kit	2250,-

LEXSY Expression Vectors

Product	Cat. No.	Amount	Price EUR
pLEXSY-ble2 Constitutive expression vector, LEXSY Bleo selection	EGE-231	5 µg	480,-
pLEXSY-hyg2 Constitutive expression vector, LEXSY Hygro selection	EGE-232	5 µg	480,-
pLEXSY-neo2 Constitutive expression vector, LEXSY Neo selection	EGE-233	5 µg	480,-
pLEXSY-sat2 Constitutive expression vector, LEXSY NTC selection	EGE-234	5 µg	480,-
pLEXSY_I-blecherry3 Inducible int. expression vector, LEXSY Bleo selection	EGE-243	5 µg	480,-
pLEXSY_I-ble3 Inducible int. expression vector, LEXSY Bleo selection	EGE-244	5 µg	480,-
pLEXSY_I-neo3 Inducible int. expression vector, LEXSY Neo selection	EGE-245	5 µg	480,-
pLEXSY_IE-blecherry4 Inducible ep. expression vector, LEXSY Bleo selection	EGE-251	5 µg	480,-

LEXSY Cultivation Kits

Product	Cat. No.	Amount	Price EUR
LEXSY Cultivation Kit P10 contains LEXSY host P10, for constitutive expression	LT-102	1 Kit	720,-
LEXSY Cultivation Kit T7-TR contains LEXSY host T7-TR, for inducible expression	LT-111	1 Kit	1650,-

LEXSY Strains

Product	Cat. No.	Amount	Price EUR
LEXSY host P10 for constitutive expression	LT-101	3x 1,6 ml	480,-
LEXSY host T7-TR for inducible expression	LT-110	3x 1.6 ml	1400,-

LEXSY Plating Kits

Product	Cat. No.	Amount	Price EUR
LEXSY Plating Kit comfort components for solid medium, LEXSY BHI- and fetal-calf-serum-based, with nitrocellulose membranes, spatula, dishes & serological pipettes	ML-451	for 40 plates	800,-
LEXSY Plating Kit core components for solid medium, LEXSY BHI- and fetal-calf-serum-based, without nitrocellulose membranes, spatula, dishes & serological pipettes	ML-452	for 40 plates	650,-
LEXSY Plating Kit basic components for solid medium, LEXSY BHI -based, without fetal calf-serum, nitrocellulose membranes, spatula, dishes & serological pipettes	ML-453	for 40 plates	630,-

Complex Media

Product	Cat. No.	Amount	Price EUR
LEXSY BHI - Liquid Media Kit sterile, brain-heart-infusion based medium, recommended for strain maintenance, electroporation, expression evaluation and cryoconservation	ML-411S	1 L	90,-
	ML-411L	5 L	360,-
LEXSY BHI - Powder Media Kit brain-heart-infusion based medium, recommended for strain maintenance, electroporation, expression evaluation and cryoconservation	ML-412S	for 1 L	45,-
	ML-412L	for 5 L	180,-
	ML-412XL	for 10 L	350,-
	ML-412XXL	for 50 L	1400,-
LEXSY YS - Liquid Media Kit sterile, yeast-soybean based, casein-free medium free of animal components	ML-431S	1 L	100,-
	ML-431L	5 L	400,-
LEXSY YS - Powder Media Kit yeast-soybean based, casein-free medium free of animal components	ML-432S	for 1 L	50,-
	ML-432L	for 5 L	200,-
	ML-432XL	for 10 L	375,-
	ML-432XXL	for 50 L	1500,-

Synthetic Media

Product	Cat. No.	Amount	Price EUR
Synthetic LEXSY Medium - liquid, ready-to-grow sterile, contains Hemin and Pen-Strep, shelf life 2 weeks	ML-103S	1 L	150,-
	ML-103L	5 L	600,-
Synthetic LEXSY - Liquid Media Kit sterile, with Hemin and Pen-Strep stock solutions, shelf life 6 months	ML-107S	1 L	150,-
	ML-107L	5 L	600,-

Additives

Product	Cat. No.	Amount	Price EUR
Hemin (porcine) sterile 500x stock solution in 50% triethanolamine (for 1, 5, 10 or 50 L of medium)	ML-108S	2 ml	10,-
	ML-108L	10 ml	40,-
	ML-108XL	20 ml	75,-
	ML-108XXL	100 ml	300,-
Pen-Strep sterile 200x stock solution of penicillin and streptomycin (for 1, 5, 10 or 50 L of medium)	ML-105S	5 ml	10,-
	ML-105L	25 ml	40,-
	ML-105XL	50 ml	75,-
	ML-105XXL	250 ml	300,-

Antibiotics

Product	Cat. No.	Amount	Price EUR
LEXSY NTC 1.000x sterile stock solution, 100 mg/ml	AB-101S	1 ml	40,-
	AB-101L	5 ml	160,-
	AB-101-10ML	10 ml	300,-
	AB-101-50ML	50 ml	1200,-
LEXSY NTC powder (non-sterile)	AB-102L	1 g	300,-
	AB-102XL	5 g	1200,-
	AB-102-25G	25 g	3650,-
LEXSY Bleo 1.000x sterile stock solution, 100 mg/ml	AB103S	1 ml	80,-
	AB-103L	5 ml	320,-
LEXSY Hygro 1.000x sterile stock solution, 100 mg/ml	AB-104S	1 ml	80,-
	AB-104L	5 ml	320,-
LEXSY Neo 1.000x sterile stock solution, 50 mg/ml	AB-105S	1 ml	40,-
	AB-105L	5 ml	160,-
LEXSY Tet 1.000x sterile stock solution, 10 mg/ml, for inducible protein expression in LEXSY host T7-TR	AB-106S	1 ml	5,-
	AB-106L	5 ml	20,-
	AB-106XL	10 ml	40,-
	AB-106XXL	50 ml	150,-