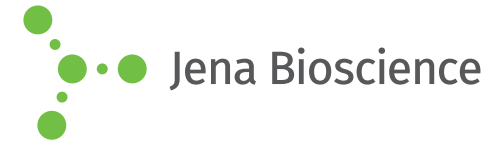




JBScreen Detergents 3

Cat.-No.: CS-523

SCREEN FORMULATION



No.	Detergent	Type	CAS Number	MW	CMC (mM)
E1	40.00 mM LysoFos Choline 10	Synthetic Lipid	22248-63-1	411.4	8.00
E2	250 mM MEGA-9	Non ionic	85261-19-4	335.4	25.00
E3	0.9 mM BRIJ® 35	Non ionic	9002-92-0	1198.0	0.09
E4	8.1 mM C10E5	Non ionic	23244-49-7	378.5	0.81
E5	40.00 mM CHAPSO	Zwitterionic	82473-24-3	630.9	8.00
E6	38.00 mM CYMAL®-4	Non ionic	181135-57-9	480.5	7.60
E7	10.8 mM CY-TRIPGLU	Non ionic	1126527-71-6	665.8	1.80
E8	18.00 mM n-Decyl-β-D-maltoside	Non ionic	82494-09-5	482.6	1.80
E9	7.2 mM Dimethyldodecylphosphine oxide	Non ionic		232.3	1.20
E10	106.4 mM FOS-Choline®-ISO-11	Zwitterionic	869647-65-4	337.4	26.60
E11	35.00 mM HEGA-10	Non ionic	139361-84-5	379.5	7.00
E12	1.8 mM Façade-EM	Non ionic		1115.3	0.18
F1	9.6 mM LAPAO	Zwitterionic	61792-31-2	300.6	1.60
F2	76.00 mM n-Decyl-N,N-dimethylglycine	Zwitterionic	2644-45-3	243.4	19.00
F3	9.00 mM n-Dodecyl-N,N-dimethylglycine	Zwitterionic	683-10-3	271.4	1.50
F4	0.3 mM n-Tridecyl-β-D-maltoside	Non ionic	93911-12-7	524.6	0.03
F5	5.9 mM n-Undecyl-β-D-maltoside	Non ionic	253678-67-0	496.6	0.59
F6	21.6 mM Ph-TRIPGLU	Non ionic	1126527-67-0	659.8	3.60
F7	40.00 mM Sodium deoxycholate	Ionic	302-95-4	414.6	4.00
F8	400 mM ZWITTERGENT® 3-10	Zwitterionic	15163-36-7	307.6	40.00
F9	40.00 mM C8E4	Non ionic	19327-39-0	306.5	8.00
F10	50.00 mM C8E6	Non ionic	4440-54-4	394.5	10.00
F11	2.00 mM ANAPOE®-X-114	Non ionic	9036-19-5	426.6	0.20
F12	6.5 mM ANAPOE®-X-305	Non ionic	9002-93-1	1526.0	0.65

