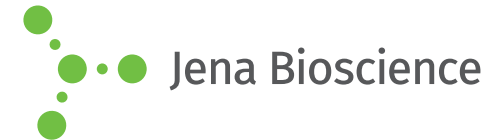




JBScreen Detergents 4

Cat.-No.: CS-524

SCREEN FORMULATION



No.	Detergent	Type	CAS Number	MW	CMC (mM)
G1	120 mM C6E4	Non ionic	39619-69-7	278.4	30.00
G2	210 mM C7E5	Non ionic	65316-79-2	336.5	21.00
G3	10.2 mM Octyl Glucose Neopentyl Glycol	Non ionic	1257853-32-9	568.7	1.02
G4	0.8 mM C12E9	Non ionic	3055-99-0	582.8	0.08
G5	2.00 mM ANAPOE®-C12E10	Non ionic	6540-99-4	627.0	0.20
G6	31.00 mM FOS-Choline®-UNSAT-11-10	Zwitterionic	121045-77-0	335.4	6.20
G7	1.9 mM CYMAL®-7	Non ionic	349477-49-2	522.5	0.19
G8	790 mM MEGA-8	Non ionic	85316-98-9	321.4	79.00
G9	13.2 mM n-decyl-β-D-glucoside	Non ionic	58846-77-8	320.4	2.20
G10	20.00 mM n-Dodecyl-N,N-dimethylamine-N-oxide	Zwitterionic	1643-20-5	229.4	2.00
G11	1.5 mM n-Dodecyl-α-D-maltoside	Non ionic	116183-64-3	510.6	0.15
G12	30.00 mM n-Nonyl-β-D-maltoside	Non ionic	106402-05-5	468.5	6.00
H1	42.5 mM n-Octyl-β-D-thiomaltoside	Non ionic	148616-91-5	470.6	8.50
H2	0.5 mM n-Undecyl-β-D-thiomaltoside	Non ionic	148565-57-5	512.7	0.05
H3	96.8 mM OHES	Non ionic	7305-30-8	206.4	24.20
H4	60.00 mM 1,2-Dihexanoyl-sn-Glycero-3-Phosphocholine	Zwitterionic	34506-67-7	453.5	15.00
H5	7.00 mM LysoFos Choline 12	Synthetic Lipid	20559-18-6	439.5	0.70
H6	96.8 mM RDHPOS	Non ionic	142741-63-7	236.4	24.20
H7	115 mM Sodium cholate hydrate	Ionic	206986-87-0	430.6	11.50
H8	82.00 mM Sodium dodecyl sulfate	Ionic	151-21-3	288.4	8.20
H9	40.00 mM ZWITTERGENT® 3-12	Zwitterionic	14933-08-5	335.6	4.00
H10	1.3 mM n-Dodecyl β-D-glucoside	Non ionic	59122-55-3	348.5	0.13
H11	2.3 mM ANAPOE® X-100	Non ionic	9002-93-1	647.0	0.23
H12	0.36 mM PCC-α-maltoside	Non ionic		547.7	0.04

