



JBScreen Detergents 1

Cat.-No.: CS-521

SCREEN FORMULATION



No.	Detergent	Type	CAS Number	MW	CMC (mM)
A1	6.00 mM Octyl maltoside, fluorinated	Non ionic	118680-70-9	688.4	1.02
A2	3.00 mM ANAPOE® NID-P40	Non ionic	2497-59-8	603.0	0.30
A3	8.4 mM 1,2-Diheptanoyl-sn-Glycero-3-Phosphocholine	Zwitterionic	39036-04-9	481.6	1.40
A4	8.4 mM Big CHAP, deoxy	Non ionic	86303-23-3	862.1	1.40
A5	0.9 mM C12E8	Non ionic	3055-98-9	538.8	0.09
A6	7.8 mM ANAPOE®-C10E9	Non ionic	26183-52-81	555.0	1.30
A7	56.00 mM CYCLOFOS®-4	Zwitterionic	675126-15-5	320.9	14.00
A8	112 mM CYGLU-3	Non ionic	869541-00-4	304.4	28.00
A9	5.6 mM CYMAL®-6	Non ionic	228579-27-9	508.5	0.56
A10	1.7 mM n-Dodecyl-β-D-maltoside	Non ionic	69227-93-6	510.6	0.17
A11	5.7 mM Dimethyldodecylphosphine oxide	Non ionic	871-95-4	246.4	0.57
A12	44.00 mM FOS-Choline®-10	Zwitterionic	70504-28-8	323.4	11.00
B1	78.00 mM HECAMEG	Non ionic	115457-93-5	335.4	19.50
B2	1250 mM n-Hexyl-β-D-glucoside	Non ionic	59080-45-4	264.4	250.00
B3	70.00 mM MEGA-10	Non ionic	85261-20-7	349.5	7.00
B4	9.6 mM n-Decyl-α-D-maltoside	Non ionic	168037-12-5	482.6	1.60
B5	25.8 mM DDMAB	Zwitterionic	15163-30-1	299.5	4.30
B6	32.5 mM n-Nonyl-β-D-glucoside	Non ionic	69984-73-2	306.4	6.50
B7	72.00 mM n-Octyl-β-D-glucoside	Non ionic	29836-26-8	292.4	18.00
B8	40.00 mM CHAPS	Zwitterionic	75621-03-3	614.9	8.00
B9	19.2 mM n-Undecyl-N,N-Dimethylamine-Oxide	Zwitterionic	15178-71-9	215.4	3.20
B10	0.12 mM TWEEN® 80	Non ionic	9005-65-6	1309.7	0.01
B11	15.00 mM n-Decanoylsucrose	Non ionic	31835-06-0	496.6	2.50
B12	1650 mM ZWITTERGENT® 3-08	Zwitterionic	15178-76-4	279.4	330.00

