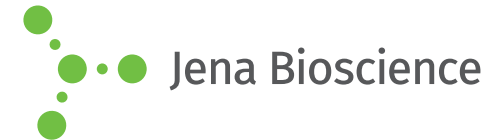




JBScreen Detergents 2

Cat.-No.: CS-522

SCREEN FORMULATION



No.	Detergent	Type	CAS Number	MW	CMC (mM)
C1	2.7 mM 1,2-dioctanoyl-sn-glycero-3-phosphocholine	Zwitterionic	19191-91-4	509.6	0.27
C2	110 mM 2,6-Dimethyl-4-heptyl-β-D-maltoside	Non ionic	869638-31-3	468.5	27.50
C3	1000 mM C6E3	Non ionic	25961-89-1	234.3	100.00
C4	35.5 mM C8E5	Non ionic	19327-40-3	350.5	7.10
C5	1.00 mM ANAPOE®-C13E8	Non ionic	9043-30-5	553.0	0.10
C6	16.08 mM CYCLOFOS®-6	Zwitterionic	657393-65-2	349.2	2.68
C7	120 mM CYMAL®-3	Non ionic	181135-58-0	466.5	30.00
C8	0.36 mM Decyl Maltose Neopentyl Glycol	Non ionic	1257852-99-5	949.8	0.04
C9	0.1 mM Lauryl Maltose Neopentyl Glycol	Non ionic	1257852-96-2	1005.2	0.01
C10	9.00 mM FOS-Choline®-12	Zwitterionic	29557-51-5	351.5	1.50
C11	30.00 mM Genapol® X-100	Non ionic	9043-30-5	641.0	7.50
C12	790 mM n-Heptyl-β-D-glucopyranoside	Non ionic	78617-12-6	278.4	79.00
D1	40.00 mM Dimethylnonylphosphine oxide	Non ionic		204.1	10.00
D2	42.00 mM n-Decyl-N,N-Dimethylamine-Oxide	Non ionic	2605-79-0	201.4	10.50
D3	9.00 mM n-Decyl-β-D-thiomaltoside	Non ionic	14565-56-4	498.6	0.90
D4	116 mM n-Heptyl-β-D-thioglucoiside	Non ionic	85618-20-8	294.4	29.00
D5	78.00 mM n-Octyl-β-D-maltoside	Non ionic	82494-08-4	454.4	19.50
D6	2.00 mM Nonidet P40	Non ionic	127087-87-0	573.0	0.20
D7	1.2 mM FOS-Choline®-14	Zwitterionic	77733-28-9	379.5	0.12
D8	46.00 mM C-HEGA-11	Non ionic	864434-16-2	391.5	3.00
D9	0.6 mM ZWITTERGENT 3-16	Zwitterionic	2281-11-0	391.7	0.06
D10	27.00 mM TRIPAO	Zwitterionic	318293-46-8	362.5	4.50
D11	0.59 mM TWEEN 20	Non ionic	9005-64-5	1228.0	0.06
D12	7.2 mM Omega-Undecylenyl-β-D-Maltoside	Non ionic	869704-76-7	494.6	1.20

