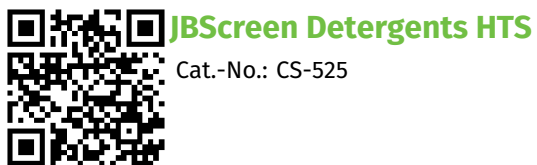


No.	Detergent	Type	CAS Number	MW	CMC (mM)
A1	6.00 mM Octyl maltoside, fluorinated	Non ionic	118680-70-9	688.4	1.02
A2	3.00 mM ANAPOE® NID-P40	Non ionic	2497-59-8	603.0	0.30
A3	8.4 mM 1,2-Diheptanoyl-sn-Glycero-3-Phosphocholine	Zwitterionic	39036-04-9	481.6	1.40
A4	8.4 mM Big CHAP, deoxy	Non ionic	86303-23-3	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

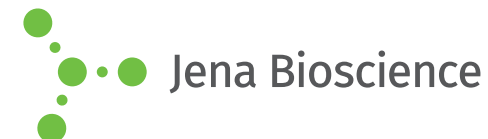
No.	Detergent	Type	CAS Number	MW	CMC (mM)
C1	2.7 mM 1,2-dioctanoyl-sn-glycero-3-phosphocholine	Zwitterionic	19191-91-4	509.6	0.27
C2	110 mM 2,6-Dimethyl-4-heptyl-β-D-maltoside	Non ionic	869638-31-3	468.5	27.50
C3	1000 mM C6E3	Non ionic	25961-89-1	234.3	100.00
C4	35.5 mM C8E5	Non ionic	19327-40-3	350.5	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

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 Dimethylundecylphosphine 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



Cat.-No.: CS-525

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

