

biological detergents

Name	Description	Brand	Cat. No.	CAS	MW (g/mol)	CMC	Aggregation Number	HLB	Cloud Point (°C)	Average Micellar Weight	Diagnostic Applications	Molecular Biology	Cell Culture	Electrophoresis/ Chromatography	Membrane Protein Solubilization	Enzymology	Antigen/Vaccine Preparation	Drug Delivery/ Liposomes	Mass Spectrometry	Microbiology
Anionic Detergents																				
1-Octanesulfonic acid sodium salt	BioXtra	Sigma-Aldrich	00133	5324-84-5	216.27															
Chenodeoxycholic acid	≥97% (TLC) from ox or sheep bile, ≥98% (TLC)	Sigma	C9377	474-25-9	392.57	3 mM														
Deoxycholic acid	BioXtra, ≥99% (HPLC) ≥ 98% (HPLC)	Sigma	C1129	81-25-4	408.57	2-6 mM	3-12	16	1,200-5,000											
Docusate sodium salt	BioXtra, ≥99.0% (TLC) BioXtra, ≥99%	Sigma	30960		444.56															
LDS, Lithium dodecyl sulfate	BioReagent, for molecular biology, suitable for electrophoresis, ≥ 98.5% (GC) BioXtra, ≥ 98.5% (GC)	Sigma-Aldrich	C1781	2044-56-6	272.33	8.2 mM	62	>100	17,000											
N-Lauroylsarcosine	BioReagent, for molecular biology, ≥ 97.0% (GC) BioXtra, ≥ 95% (titration)	Sigma	14632		271.4	14.6 mM	2		600											
N-Lauroylsarcosine sodium salt	purum p.a., ≥ 98.0% (GC) BioUltra, for molecular biology, ≥ 99.0% (HPLC) BioXtra, ≥ 97% (TLC) ≥ 98% (titration)	Sigma	61747	137-16-6	293.38	14.6 mM	2		600											
Sodium 1-decanesulfonate	BioReagent, for molecular biology, ≥94% 20% solution, for molecular biology	Sigma	17414		244.33															
Sodium cholate hydrate	BioReagent, suitable for cell culture, ≥99% (HPLC) ≥97.0% (dried material, NT) BioXtra, ≥99% (HPLC) from ox and/or sheep bile, ≥99% (HPLC)	Sigma	27029		430.55	9-15 mM	2-3	18	900-1,300											
Sodium deoxycholate	BioXtra, ≥98.0% (dry matter, NT) ≥97% (titration)	Sigma	30970	302-95-4	414.55	2-6 mM	3-12	16	1,200-5,000											
Sodium deoxycholate monohydrate	BioUltra, for manufacturing of diagnostic kits and reagents BioXtra, ≥99.0% (NT)	Sigma	546046		432.57	2-6 mM	3-12	16	1,200-5,000											
SDS, Sodium dodecyl sulfate	BioUltra, for molecular biology, 10% in H ₂ O ≥ 98.0% (GC) BioUltra, for molecular biology, ≥ 99.0% (GC) BioUltra, for molecular biology, 20% in H ₂ O BioReagent, for molecular biology, suitable for electrophoresis, ≥ 98.5% (GC) ≥ 90% ACS reagent, ≥ 99.0% (titration) 92.5-100.5% based on total alkyl sulfate content ReagentPlus™, ≥ 98.5% (GC) BioXtra, ≥ 99.0% (GC) ≥99.0% (GC), dust-free pellets Testing conforms to Pharmacopoeia anhydrous, free-flowing, Redi-Dri™, ACS reagent, ≥99.0%	Sigma	30033	13419-61-9	244.33															
Sodium glycocholate	≥97% (HPLC)	Sigma	60789	16564-43-5	471.61															
Sodium glycocholate hydrate	≥97% (TLC)	Sigma	C7132	138950-81-5	487.6	13 mM	1.9		900											
Sodium taurocholate	BioXtra, ≥99% (TLC)	Sigma	C0910	16409-34-0	471.61	2.1 mM	2		900											
Sodium taurodesoxycholate	≥98% (elemental analysis)	Sigma-Aldrich	15249	2832-45-3	188.22															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Cationic Detergents																				
Alkyltrimethylammonium bromide	≥95% (TLC)	Sigma	M7635	8044-71-1																
Benzalkonium chloride	BioXtra	Sigma	86295	63449-41-2	396.09															
Benzyltrimethylammonium chloride	cationic detergent	Sigma	94136	122118-0																
Dimethyldioctadecylammonium bromide	≥98% (TLC)	Sigma	D2779	3700-67-2	630.95															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Akytrimethylammonium bromide	≥95% (TLC)	Sigma	M7635	8044-71-1																
Benzalkonium chloride	BioXtra	Sigma	86295	63449-41-2	396.09															
Benzyltrimethylammonium chloride	cationic detergent	Sigma	94136	122118-0																
Dimethyldioctadecylammonium bromide	≥98% (TLC)	Sigma	D2779	3700-67-2	630.95															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Akytrimethylammonium bromide	≥95% (TLC)	Sigma	M7635	8044-71-1																
Benzalkonium chloride	BioXtra	Sigma	86295	63449-41-2	396.09															
Benzyltrimethylammonium chloride	cationic detergent	Sigma	94136	122118-0																
Dimethyldioctadecylammonium bromide	≥98% (TLC)	Sigma	D2779	3700-67-2	630.95															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Akytrimethylammonium bromide	≥95% (TLC)	Sigma	M7635	8044-71-1																
Benzalkonium chloride	BioXtra	Sigma	86295	63449-41-2	396.09															
Benzyltrimethylammonium chloride	cationic detergent	Sigma	94136	122118-0																
Dimethyldioctadecylammonium bromide	≥98% (TLC)	Sigma	D2779	3700-67-2	630.95															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Akytrimethylammonium bromide	≥95% (TLC)	Sigma	M7635	8044-71-1																
Benzalkonium chloride	BioXtra	Sigma	86295	63449-41-2	396.09															
Benzyltrimethylammonium chloride	cationic detergent	Sigma	94136	122118-0																
Dimethyldioctadecylammonium bromide	≥98% (TLC)	Sigma	D2779	3700-67-2	630.95															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Akytrimethylammonium bromide	≥95% (TLC)	Sigma	M7635	8044-71-1																
Benzalkonium chloride	BioXtra	Sigma	86295	63449-41-2	396.09															
Benzyltrimethylammonium chloride	cationic detergent	Sigma	94136	122118-0																
Dimethyldioctadecylammonium bromide	≥98% (TLC)	Sigma	D2779	3700-67-2	630.95															
Sodium pentanesulfonate	BioXtra (elemental analysis)	Sigma-Aldrich	H9026	22767-49-3	174.19															
Sodium taurocholate hydrate	BioXtra, ≥95% (TLC) ≥ 97.0% (TLC) ≥ 95% (TLC)	Sigma	19034	345909-26-4	537.68	3-11 mM	4		2,100											
Sodium taurodesoxycholate hydrate	BioXtra, ≥ 97% (TLC)	Sigma	10557	207737-97-1	521.69	1-4 mM	6		3,100											
Sodium tauroursodeoxycholate	≥ 95% (TLC)	Sigma	10266	14605-22-2	521.69	4 mM		20.1												
Non-Ionic Detergents																				
1-Oleoyl- α -D-glycerol	≥99.0% (TLC)	Sigma	M7765	1111-03-5	356.54															
Brij® L23	Solution, 30 % (w/v) in H ₂ O suitable for Stern-Moore chromatography main component: tricosaethylene glycol dodecyl ether	Sigma-Aldrich	84184	9002-92-0	1198.0 (avg.)	0.091 Mm	40	16.9	>100	48,000										
Brij® O10	non-ionic surfactant	Sigma	86136	9004-98-2	709.00				12.4											