



Promega

Special Offer

SALE!

Great Offers on
High Quality
Biochemicals
Products

**WE'RE HERE
FOR YOU**



SAVE

NOW

Offer Valid From 10th Feb to March 31st 2023

Promotion Code **PROMEGA_FM2023**



Save 20 %**BioChemicals**

Part Number	Product Name	Size	INR Price 2022-23	Offer Price
D1501	Lambda DNA	250microg	7480	5984
D1811	Herring Sperm DNA	10mg	4845	3876
D1815	Herring Sperm DNA	100mg	13685	10948
G1881	Blue/Orange Loading Dye, 6X	3ml	3740	2992
H1181	Diamond Nucleic Acid Dye	500microl	7055	5644
H5003	Boric Acid, Molecular Biology Grade	1kg	7480	5984
H5032	EDTA, Disodium Salt, Molecular Biology Grade	500g	9945	7956
H5041	Ethidium Bromide Solution, Molecular Grade	10ml	3995	3196
H5052	Formamide, Molecular Grade	500 ml	10965	8772
H5073	Glycine, Molecular Biology Grade	1kg	9945	7956
H5114	Sodium Dodecyl Sulfate, Molecular Biology Grade (SDS)	500 g	13685	10948
H5113	Sodium Dodecyl Sulfate, Molecular Biology Grade (SDS)	100g	5100	4080
H5123	Tris-HCl, Molecular Biology Grade	500g	10115	8092
H5125	Tris-HCl, Molecular Biology Grade	2,500g	41140	32912
H5131	Tris Base, Molecular Biology Grade	500g	6545	5236
H5135	Tris Base, Molecular Biology Grade	2,500g	23460	18768
H5141	Triton X-100, Molecular Biology Grade	500ml	9775	7820
H5151	Tween 20, Molecular Biology Grade	500ml	7565	6052
H5273	Sodium Chloride, Molecular Biology Grade	1kg	5100	4080
H5302	HEPES, Molecular Biology Grade (free acid)	100g	7140	5712
H5383	Guanidine-HCl, Molecular Biology Grade	500g	15045	12036
H5433	Glycerol, Molecular Biology Grade	1,000ml	5270	4216
P1193	Nuclease-Free Water	50ml	4428	3542
V2111	Agarose, Low Melting Point, Analytical Grade	25g	13430	10744
V2791	Guanidine Thiocyanate, Molecular Grade	100g	4505	3604
V3011	PEG 8000 Powder, Molecular Biology Grade	500g	6715	5372
V3021	Proteinase K	100mg	14500	11600
V3115	Acrylamide, Molecular Grade	500g	14025	11220
V3143	Bisacrylamide, Molecular Grade	125g	9010	7208
V3151	DTT, Molecular Grade (Dry Powder)	5g	7905	6324
V3155	DTT, Molecular Grade (Dry Powder)	25g	28475	22780
V3171	Urea	1kg	5865	4692
V3175	Urea	5kg	19295	15436
V3841	Agarose, LMP, Preparative Grade for Small Fragments (10 to 1,000bp)	25g	12665	10132
V3941	X-Gal	100mg/2ml	7650	6120
V3951	IPTG, Dioxane-Free	5g	18275	14620
V3953	IPTG, Dioxane-Free	50g	107780	86224
V7983	Antibiotic G-418 Sulfate	5g	41820	33456

Terms & Conditions

Offer is valid from 10th Feb, 2023 till 31st March, 2023. Promotion Code PROMEGA_FM2023 when placing your order. Promega Biotech India Private Limited reserves the right to change or cancel the offer at anytime without any prior notice. Valid only on the products listed and cannot be combined with any other special offer, promotion, or price discount. Not valid on any prior orders. Not valid with any other offer, contract or agreement. Void where prohibited by law. Rate contract discount not applicable. Prices in Indian Rupees and is available only in India. Taxes Extra as Applicable.

Contact US :-**Promega Biotech India Private Limited**

Unit No S -114, Salcon Aurum Building Jasola District Centre, New Delhi-110044

Tel +91 11 43005817 Email : ind_custserv@promega.com

