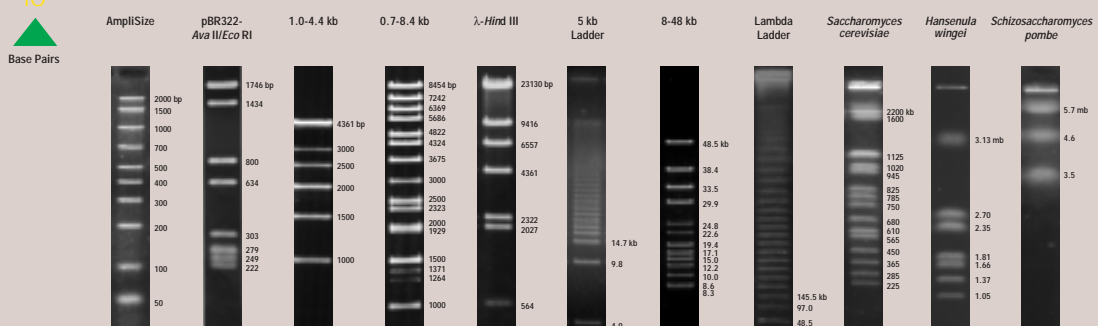
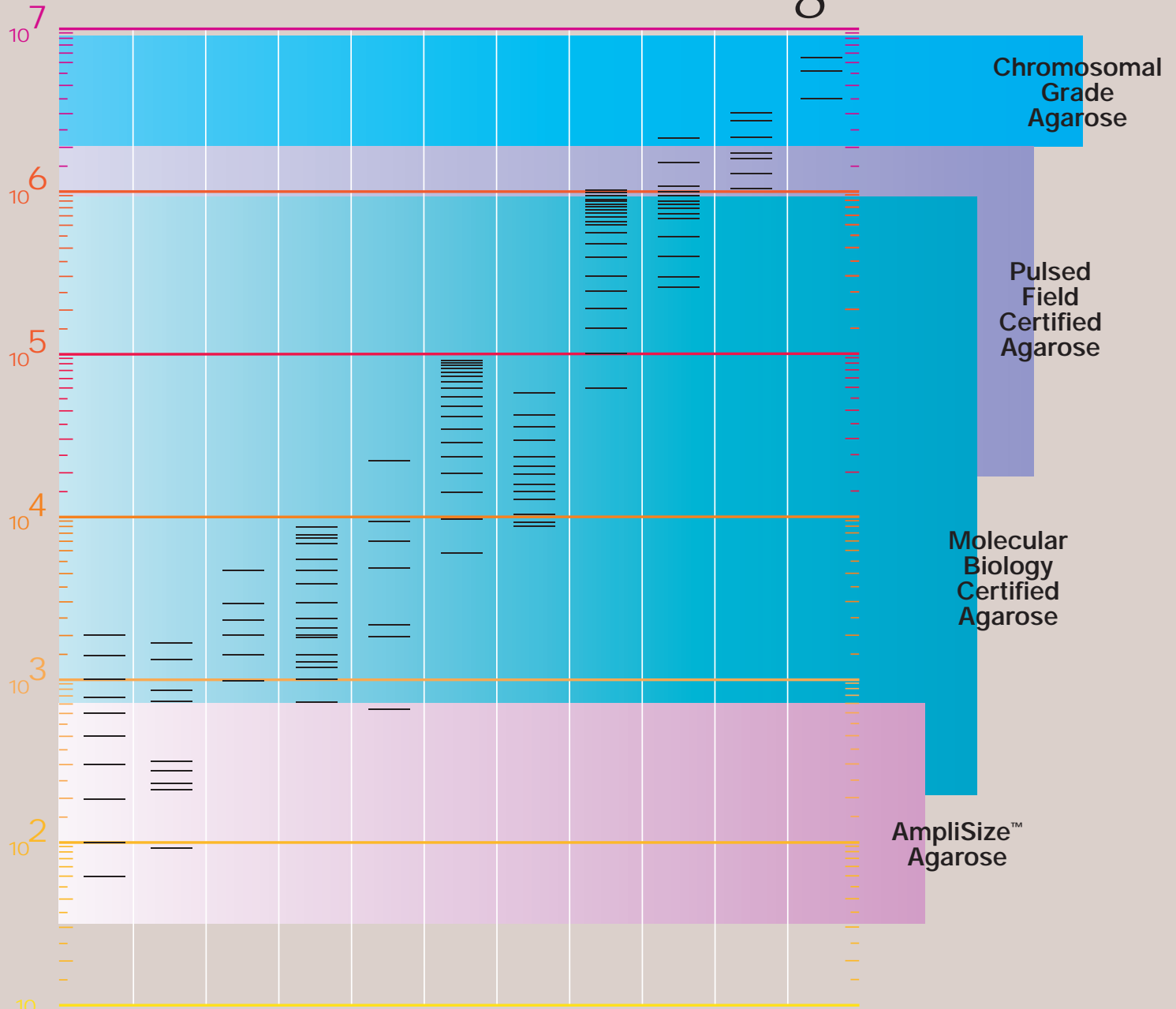




DNA Size Standards and Agaroses



Ordering Information			
Catalog No.	Product Description	Amount	Price
170-8200	Amplisize™	50 µg/pt	95.00
170-3465	pBR322-Ava II/Eco RI	10 µg	82.00
170-8210	1.0-4.4 kb ladder	50 µg/pt	95.00
170-8220	0.7-8.4 kb	100 µg	95.00
170-3470	λ-Hind III	200 µg	82.00
170-3424	5 kb Ladder	20 µg	98.00
170-3707	8-48 kb	25 µg	50.00
170-3435	λ Ladder	5 blocks	82.00
170-3605	Saccharomyces cerevisiae	5 blocks	82.00
170-3667	Hansenula wingei	5 blocks	82.00
170-3633	Schizosaccharomyces pombe	5 blocks	82.00
162-0144	Amplisize™	50 g	149.00
162-0133	Molecular Biology Cert.	100 g	115.00
162-0134		500 g	465.00
162-0137	Pulsed Field Certified	100 g	99.00
162-0138		500 g	445.00
162-0135	Chromosomal Grade	100 g	99.00
162-0136		500 g	360.00
162-0017	Low Melt Preparative	25 g	99.00
162-0019		100 g	325.00

Amplisize Agarose: Amplisize Standard, pBR322-Ava II/Eco RI Molecular Biology Certified Agarose: 1.0-4.4 kb, 0.7-8.4 kb, λ-Hind III, 5 kb ladder Pulsed Field Certified Agarose: 8-48 kb, lambda ladder, S. cerevisiae Chromosomal Grade Agarose: H. wingei, S. pombe