

Ordering Information

<i>Cat. No.</i>	<i>Product Description</i>
170-8201	20 bp Molecular Ruler DNA Size Standard
170-8202	100 bp Molecular Ruler DNA Size Standard
170-8206	100 bp PCR Ruler DNA Size Standard
170-8203	500 bp Molecular Ruler DNA Size Standard
170-8204	1 kb Molecular Ruler DNA Size Standard
170-8205	2.5 kb Molecular Ruler DNA Size Standard
170-8207	Mass Ruler DNA Mass Standard

**The Polymerase Chain Reaction (PCR) process is covered by patents owned by Hoffmann-LaRoche. Use of the PCR process requires a license.*

BIO-RAD

**Bio-Rad
Laboratories**

Molecular Bioscience Group

U.S. (800) 4BIORAD **California** (510) 741-1000 **Australia** 02-9914-2800 **Austria** (1) 877 89 01 **Belgium** 09-385 55 11 **Canada** (905) 712-2771 **China** (86-10) 2046622
Denmark 39 17 9947 **Finland** 90 804 2200 **France** (1) 43 90 46 90 **Germany** 089 31884-0
India 91-11-461-0103 **Italy** 02-21609 1 **Japan** 03-5811-6270 **Hong Kong** 7893300
The Netherlands 031318-540666 **New Zealand** 09-443 3099 **Singapore** (65) 272-9877
Spain (91) 661 70 85 **Sweden** 46 (0) 8 627 50 00 **Switzerland** 01-809

Bulletin 2152 US/EG Rev A 96-5229 197 SIG 1797 Printed in USA

Molecular Rulers™
by Which All
Other Standards
are Measured



BIO-RAD

Bio-Rad Molecular Rulers Simplify Gel Analysis

Bio-Rad's tradition as the leader in precision electrophoresis continues with seven new DNA ladders. Molecular, PCR* and Mass Rulers now make it easier to analyze gels. Ladders are available in 20 bp, 100 bp, 500 bp, 1 kb, 2.5 kb and a special 100 bp PCR

Ruler. To precisely measure samples, most of these ladders include a visually-distinct reference band. The 2.5 kb ladder, a marker ideal for FIGE gels, has a band at 10 kb and extends up to 35 kb. The 100 bp PCR Ruler extends to 3 kb in exact 100 bp increments for more accurate sizing of longer PCR products. The Mass Ruler has five coordinated bands of decreasing mass: 1,000 bp containing 100 ng; 700 bp containing 70 ng; 500 bp containing 50 ng; 200 bp containing 20 ng; 100 bp containing 10 ng.

Specifications

Product Name	Range	Quantity	Applications	Shelf Life	Shipping
20 bp Molecular Ruler	20-1,000 bp, 50 bands; ref band at 200 bp	250 μ l	100	1 year @ 4 °C	RT
100 bp Molecular Ruler	100-1,000 bp, 10 bands	250 μ l	100	1 year @ 4 °C	RT
100 bp PCR Ruler	100-3,000 bp, 30 bands; ref bands at 1,000 and 3,000 bp	200 μ l	100	1 year @ 4 °C	RT
500 bp Molecular Ruler	500-8,000 bp, 16 bands; ref band at 5,000 bp	200 μ l	100	1 year @ 4 °C	RT
1 kb Molecular Ruler	1 kb-15 kb, 15 bands; ref band at 5 kb	200 μ l	100	1 year @ 4 °C	RT
2.5 kb Molecular Ruler	2.5 kb-35 kb, 15 bands; ref band at 10 kb	400 μ l	100	1 year @ 4 °C	RT
Mass Ruler	100 bp-1,000 bp (10 ng-100 ng)	250 μ l	100	1 year @ 4 °C	RT