

# 200 bp DNA Ladder

**Code No.** 3410A      **Size:** 250  $\mu$ l  
(for 100 lanes)  
**Conc.:** 200 ng/ $\mu$ l

**Supplied Reagents:**  
**6X Loading Buffer** 1 ml

**Form:**  
10 mM Tris-HCl, pH 8.0  
1 mM EDTA

**Storage:** -20°C  
(Once opened, 6X Loading Buffer should be stored at room temperature.)  
\* 2 years from date of receipt under proper storage conditions.

**Description:**  
This marker consists of 10 fragments between 200 bp and 2 kb in multiples of 200 bp and additional fragments at 2.5 kb, 3 kb, and 4 kb. The 1 kb and 2 kb fragments are brighter than the other fragments. They are all double-stranded DNA fragments.

**Application:**  
Used as a DNA molecular size marker in agarose gel electrophoresis. After dephosphorylation of this product, it can be radioactively labeled by the kinase phosphorylation reaction using MEGALABEL™ (Cat. #6070).\*  
\* Not available in all geographic locations. Check for availability in your area.

**6X Loading Buffer (Store at RT after used):**  
36% Glycerol  
30 mM EDTA  
0.05% Bromophenol Blue  
0.035% Xylene Cyanol

Add 1/5 volume of 6X Loading Buffer with DNA solution to apply on agarose gel electrophoresis. In case precipitates generated during the storage at room temperature, dissolve in warm bath before use.

#### Application example:

200 bp DNA Ladder	2.5 $\mu$ l (500 ng)
Sterile purified water	2.5 $\mu$ l
6X Loading Buffer	1 $\mu$ l

Run agarose gel electrophoresis using ex.  
1% Agarose L03 [TAKARA] (Cat. #5003),  
or 3% PrimeGel™ Agarose PCR-Sieve (Cat. #5810A).  
Perform staining with EtBr, or SYBR® Green I Nucleic Acid Gel Stain (Cat. #5760A/5761A).\*

\* Not available in all geographic locations. Check for availability in your area.

SYBR is a registered trademark of Life Technologies Corporation. MEGALABEL and PrimeGel are trademarks of Takara Bio Inc.

#### Note

This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc. Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from Takara Bio Inc. If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at [www.takara-bio.com](http://www.takara-bio.com). Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements. All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.

