

TEST IT, YOU WILL LOVE IT!



SMARTLADDER

MOLECULAR WEIGHT DNA MARKER

The SmartLadder is among the most popular ready-to-use molecular weight DNA marker, especially designed for easy DNA quantification as well as size determination. The SmartLadder is particularly suitable for routine use and cloning applications.

Sample reference: SmartLadder: MW-1700-SA (50 μ l)

Sample content: 50 μ l of SmartLadder (enough volume to perform 10 tests using a standard loading of 5 μ l)

Shipping conditions: Shipped at room temperature

Convinced by our SmartLadder?

Please visit www.eurogentec.com or contact your local account manager to receive a quotation and make huge savings on your next DNA ladder order.



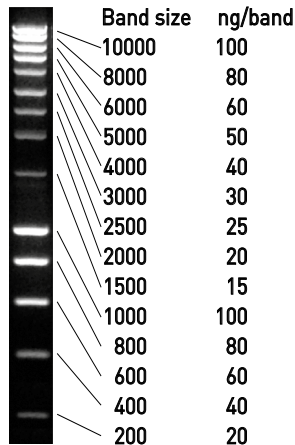
SmartLadder (200 bp – 10 kb)

The popular ready-to-use molecular weight marker for routine use

- > 1 kb ladder
- > Ready to use
- > Smart identification
- > Smart sizing
- > Smart quantitation
- > Smart storage

Sizing

14 regularly spaced bands from 200 bp to 10 kb. The 1 kb and 10 kb bands have a higher intensity than the others to allow quick and easy identification. The size of each band is a multiple of 100 bp. Each band corresponds to an exact quantity of DNA (from 15 to 100 ng).



Storage

Stable for **1 month at room temperature** and for 6 months at 4 °C. Long term storage at -20 °C.