

TEST IT, YOU WILL LOVE IT!



SMARTLADDER SF

MOLECULAR WEIGHT DNA MARKER

The SmartLadder SF is among the most popular ready-to-use molecular weight DNA marker, especially designed for easy DNA quantification as well as size determination. SmartLadder SF is perfectly suited to PCR products and short DNA fragments analysis.

Sample reference: SmartLadder SF: MW-1800-SA (50 μ l)

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

Shipping conditions: Shipped at room temperature

Convinced by our SmartLadder SF?

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 SF (100 bp - 1 kb)

The most suited molecular weight marker for PCR products analysis

- > 100 bp ladder
- > Ready to use
- > Smart identification
- > Smart sizing
- > Smart quantitation
- > Smart storage

Sizing

10 regularly spaced bands from 100 to 1000 bp. The size of each band is a multiple of 100 bp.



Storage

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

