



## CE Grade SYBR<sup>®</sup> Green I Fluorescent Dye

**Catalog Number 148-5100**

SYBR Green I is an intercalating monomeric dye with the following attractive properties: excitation maximum close to the 488 nm line of the argon ion laser, extremely low intrinsic fluorescence, and very high affinity for dsDNA.

<b>Contents</b>	1 vial CE Grade SYBR Green I Fluorescent Dye
<b>Quantity</b>	200 µl supplied in DMSO
<b>Storage</b>	Since the fluorescent dye is light sensitive, do not leave exposed to light when not in use. Store the fluorescent dye at 4 °C or below.
<b>Use</b>	Refer to the instruction manuals for the CE dsDNA Fluorescent Detection and Analysis Kits.

SYBR Green I is a registered trademark of Molecular Probe Inc., (M.P.I.). SYBR Green I is licensed from M.P.I. for CE research.