



CE Grade SYBR[®] 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.

Contents	1 vial CE Grade SYBR Green I Fluorescent Dye
Quantity	200 µl supplied in DMSO
Storage	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.
Use	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.