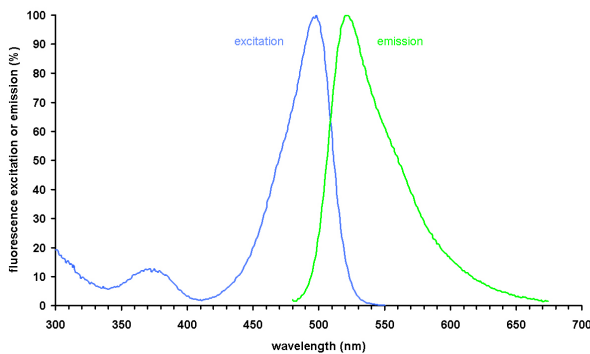




SYBR® Green Fluorescent DNA Stain

DNA intercalation dye for real-time PCR analysis

Cat. No.	Amount
PCR-378	500 µl x 100 µM
PCR-378-1ML	1 ml x 100 µM
PCR-378-10ML	10 ml x 100 µM
PCR-378-100ML	100 ml x 100 µM



Excitation (blue) and emission (green) spectrum of SYBR® Green bound to dsDNA

Description:

SYBR® Green Fluorescent DNA Stain is a superior DNA intercalator dye specially developed for DNA analysis applications including real-time PCR (qPCR). Upon binding to DNA, the non-fluorescent dye becomes highly fluorescent while showing no detectable inhibition to the PCR process. The dye is extremely stable both thermally and hydrolytically, providing convenience during routine handling.

SYBR® Green Fluorescent DNA Stain is supplied as 100 µM concentration. Vortex SYBR® Green Fluorescent DNA Stain thoroughly prior to its use. An SYBR® Green concentration of 0.5-1.0 µM in the final assay is recommended. Add SYBR® Green Fluorescent DNA Stain as indicated in the table below per assay. Please note that the preparation of a master mix may be crucial in quantitative PCR reactions to reduce pipetting errors.

Select the optical setting for SYBR® Green or FAM on the detection instrument.

final SYBR® Green concentration	20 µl PCR assay	50 µl PCR assay
0.5 µM	0.1 µl	0.25 µl
1.0 µM	0.2 µl	0.50 µl

SYBR® is a registered trademark of Invitrogen Corporation, Carlsbad, California, USA.

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: store dark

Shelf Life: 12 months

Form: liquid, supplied in 20 mM Tris-HCl pH 8.5, 0.1 mM EDTA and 0.01 % Tween-20

Color: orange

Concentration: 100 µM

Spectroscopic Properties: λ_{exc} 495 nm, λ_{em} 520 nm (bound to DNA)