

SYBR DNA Stain

Staining dye for DNA gel electrophoresis

SYBR DNA Stain is a superior DNA staining dye for DNA analysis in gel electrophoresis applications. Upon binding to DNA, the non-fluorescent dye becomes highly fluorescent allowing the detection of lowest DNA amounts. The product provides highest convenience during routine handling and avoids commonly used gel staining procedures with Ethidium Bromide.

Ordering Information

[Order Online »](#)

JBS-PCR-273	5x1.8ml
-------------	---------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance orange liquid

Emission Maximum 521 nm

Excitation Maximum 495 nm

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com