

# 5' Cy3 / 3' BHQ-2® (20 to 29 nmol)

A yield of 1 OD represents approx. 33 µg of single-stranded DNA with an equal distribution of the four bases. This corresponds to approx. 5 nmol (50 µl / 100 µM) of a DNA probe consisting of 20 bases.

## Ordering Information

[Order Online »](#)

JBS-FP-125-20	20nmol
---------------	--------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping                      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

Couple Type                Dye

Emission Maximum        563 nm

Excitation Maximum       546 nm



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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