

DEEPSEE® IGH Break Apart FISH Probe

Ready-to-use directly labeled FISH probe for
IGH Break Apart detection

The DEEPSEE® IGH Break Apart FISH Probe is a directly labeled, ready-to-use fluorescence in situ hybridization (FISH) probe for the detection of chromosomal copy number alterations and/or structural rearrangements at the IGH locus. The probe is supplied as a pre-diluted hybridization solution in formamide-based buffer and is compatible with automated hybridization instruments (HYBrite™ / ThermoBrite®). This break-apart probe flanks the IGH gene with Green / Orange labeled sequences. Structural rearrangements at this locus are indicated by separation of the two signal colors. SEEBRIGHT® Mounting Medium with DAPI is an antifade reagent that minimizes photobleaching. Use of this mounting medium significantly extends the lifetime of fluorescent signals. The inclusion of DAPI allows for staining of the nucleus. Supplied as 20 tests (200 µL) of ready-to-use probe solution.

Fluorescence in situ hybridization (FISH) is a molecular cytogenetic technique that uses fluorescently labeled DNA probes to detect and localize specific DNA sequences on chromosomes. FISH is widely used in research and clinical laboratories for the detection of chromosomal copy number changes, gene amplifications, deletions, and structural rearrangements associated with cancer and genetic disorders. DEEPSEE® directly labeled FISH probes are supplied as pre-diluted hybridization solutions, enabling high-sensitivity detection without additional labeling steps. SEEBRIGHT® Mounting Medium with DAPI is recommended for use with this probe.

Ordering Information

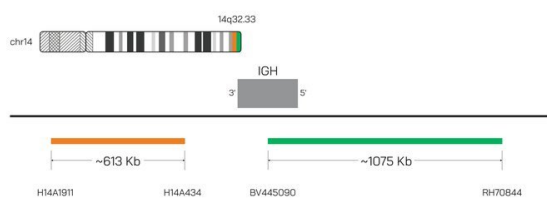
[Order Online »](#)

ENZ-GEN452-0020	20 tests
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Directly labeled, ready-to-use probe
- 2-hour hybridization; same-day results
- US manufactured with 24–48 hour delivery for uninterrupted lab workflow



Handling & Storage

Use/Stability	When used and stored as directed, reagents are stable for 1 year from date of shipment as per the date stated on the product label. Avoid freeze-thaw cycles.
Handling	Protect from light and humidity at all times. Store at -20°C. Avoid freeze-thaw cycles. Centrifuge briefly before use.
Short Term Storage	4 °C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	FISH
Assay Time	3 hr Total run time (2-hour hybridization); same-day results
Contents	1 × DEEPSEE® IGH Break Apart FISH Probe (ENZ-GEN452-0020), 20 tests (200 µL) ready-to-use probe solution in formamide-based buffer Dye Specifications: Probes labeled with Red-dUTP (Ex 580 nm/ Em 599 nm) Probes labeled with Green-dUTP (Ex 491 nm/ Em 516 nm) Probes labeled with Orange-dUTP (Ex 552 nm/ Em 576 nm) Probes labeled with Gold-dUTP (Ex 525 nm/ Em 551 nm) Probes labeled with Aqua-dUTP (Ex 434 nm/Em 480 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-
fr@enzolifesciences.com](mailto:info-fr@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)