

DEEPSEE®

PML/RARA Extended FISH Probe

Ready-to-use directly labeled FISH probe for PML/RARA Extended detection

The DEEPSEE® PML/RARA Extended 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 PML/RARA locus. The probe is supplied as a pre-diluted hybridization solution in formamide-based buffer and is compatible with automated hybridization instruments (HYBrite™ / ThermoBrite®). The probe hybridizes to the PML (Orange) and RARA (Green) loci simultaneously. Translocation events are identified by co-localization of the two signals. 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.

Ordering Information

[Order Online »](#)

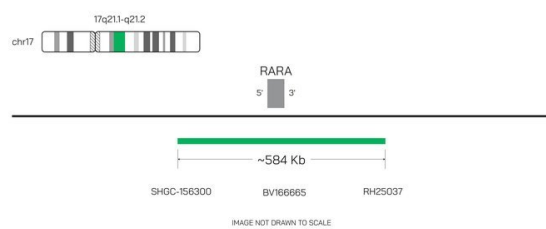
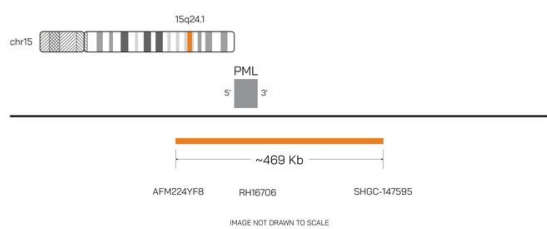
ENZ-GEN470-0005	5 tests
ENZ-GEN470-0020	20 tests

Manuals, SDS & CofA

[View Online »](#)

- Directly labeled, ready-to-use probe
- Same-day results with just 2-hour hybridization
- Uninterrupted lab workflow; 24–48 hour delivery

Compatible with both Manual and Automated (HYBrite™ / ThermoBrite®) protocols



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® PML/RARA Extended FISH Probe (ENZ-GEN470-0020), 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)