

DIGX[®] Mouse anti-digoxigenin linker (Ready-to-Use)

Ready-to-use linker for ISH applications

The DIGX[®] Mouse anti-digoxigenin linker is a ready-to-use signal amplifying reagent optimized to increase the detection sensitivity of Enzo's DIGX[®] HPV probes. Although the DIGX[®] Mouse anti-digoxigenin linker was created to enhance Enzo's DIGX[®] probes, it will recognize any digoxigenin associated nucleotide or macromolecule.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

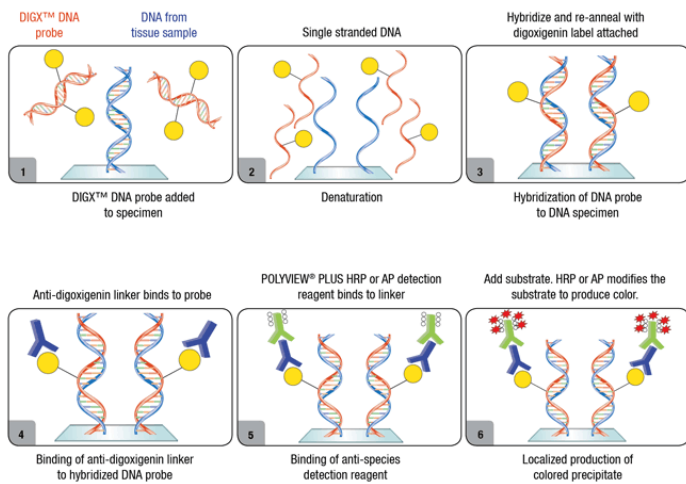
Ordering Information

[Order Online »](#)

ENZ-ABS304-6000	6ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Workflow of hybridization with DIGX[®] HPV Probes and detection with DIGXTM linkers and POLYVIEW[®] PLUS reagents

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored as recommended
Handling	Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	ISH (in situ hybridization)
Clone	HY-A.1
Formulation	Liquid. Ready-to-use in PBS containing stabilizers and preservatives.
Host	Mouse
Isotype	IgG1
Source	Purified IgG fraction of mouse serum.
Species Reactivity	Species independent
Technical Info / Product Notes	The DIGX [®] Mouse anti-digoxigenin linker can be used in conjunction with POLYVIEW [®] PLUS reagents and kits to visualize digoxigenin-labelled targets.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)