

Mouse anti-Rabbit Linker (Ready-to-use)

Ready-to-use linker for IHC applications

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

Ordering Information

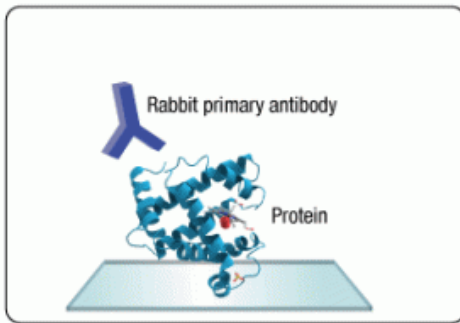
[Order Online »](#)

ENZ-ACC115-6000	6ml
-----------------	-----

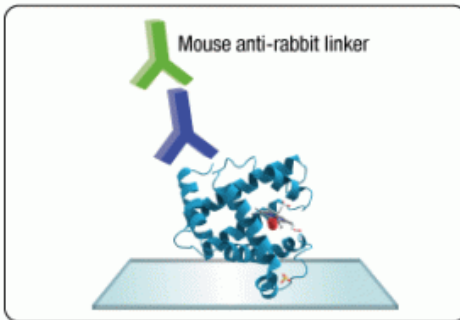
Manuals, SDS & CofA

[View Online »](#)

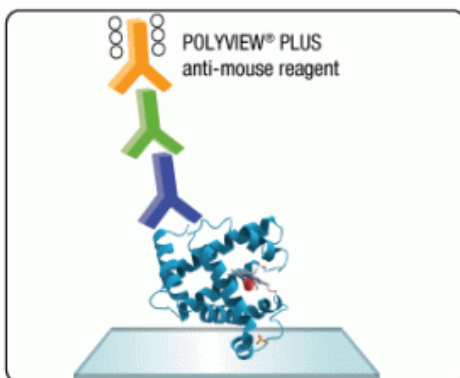
1. Add Rabbit primary antibody



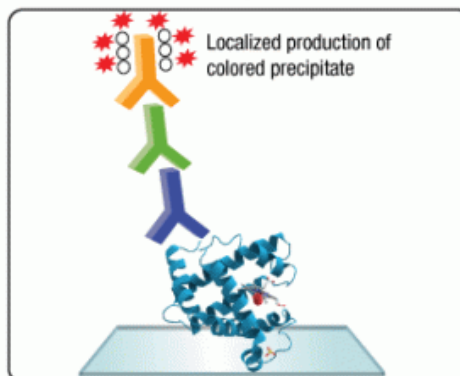
2. Add Mouse anti-rabbit linker



3. POLYVIEW® PLUS anti-mouse reagent binds to linker



4. Add Substrate HRP or AP modifies the substrate to produce color



Workflow of Mouse anti-Rabbit Linker and subsequent detection with POLYVIEW® PLUS (anti-Mouse) 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	IHC, ISH (in situ hybridization)
Application Notes	The Mouse anti-Rabbit Linker can be used in conjunction with POLYVIEW [®] PLUS (anti-Mouse) Reagents and kits to visualize rabbit targets.
Formulation	Ready-to-use solution in PBS containing stabilizers and preservatives
Host	Mouse
Isotype	IgG
Species Reactivity	Rabbit
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)