

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

Ready-to-use linker for IHC applications

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

## Ordering Information

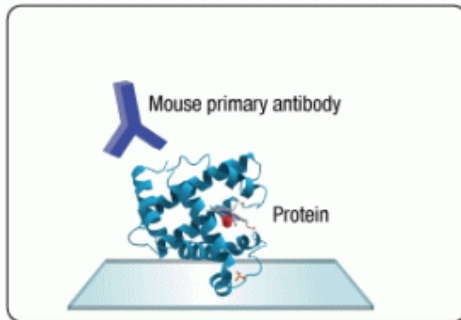
[Order Online »](#)

ENZ-ACC116-6000	6ml
-----------------	-----

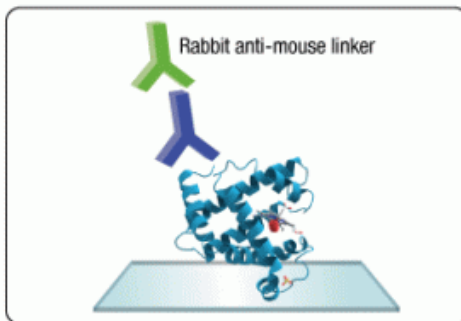
## Manuals, SDS & CofA

[View Online »](#)

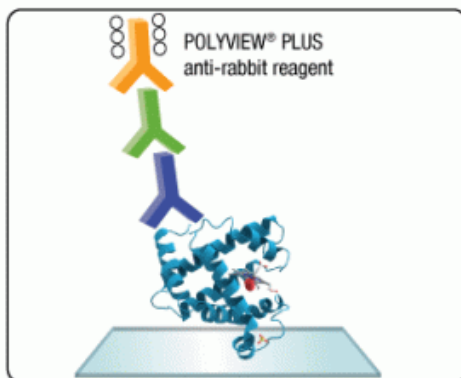
**1. Add Mouse primary antibody**



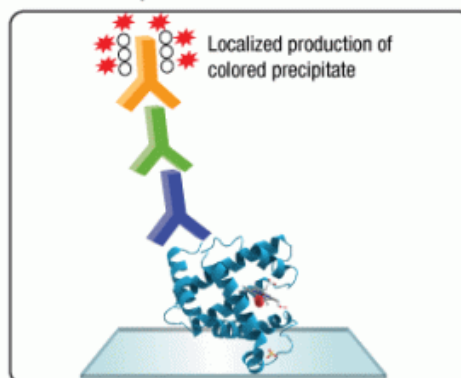
**2. Add Rabbit anti-mouse linker**



**3. POLYVIEW® PLUS anti-rabbit reagent binds to linker**



**4. Colored reaction precipitate produced with substrate**



## 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 Rabbit anti-Mouse Linker can be used in conjunction with POLYVIEW <sup>®</sup> PLUS (anti-Rabbit) Reagents and kits to visualize mouse targets.
Formulation	Ready-to-use solution in PBS containing stabilizers and preservatives
Host	Rabbit
Isotype	IgG
Species Reactivity	Mouse
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)