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 »

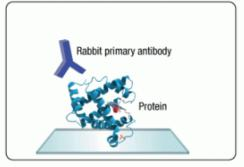
ENZ-ACC115-6000

6ml

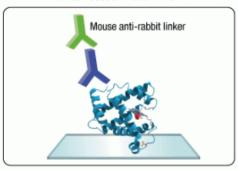
Manuals, SDS & CofA

View Online »

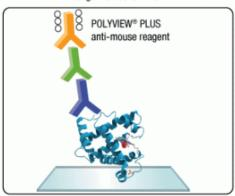




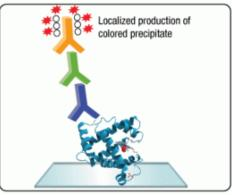
2. Add Mouse anti-rabbit linker



 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 NotesThe 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

.

Last modified: May 29, 2024

