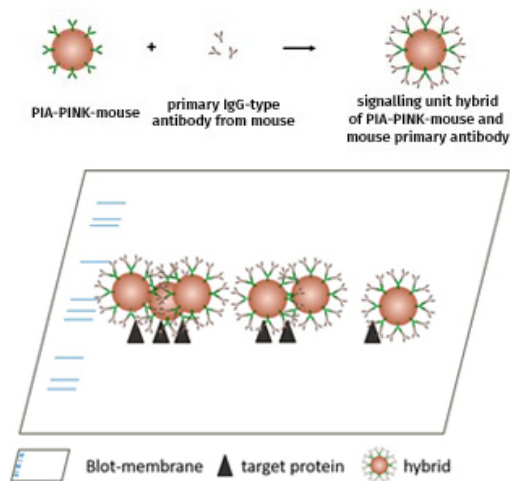




Particle Immuno Assay PIA-PINK-mouse

Particle Immuno Assay
2nd IgG antibody nanoparticles for Immunoassay

Cat. No.	Amount
PS-112	5 Blots



Description:

The PIA-PINK staining solution is a nanoparticle of gold-conjugated anti-rabbit or anti-mouse 2nd antibodies. When the antigen-specific primary antibody (protein-specific primary antibody) is added by the user, the secondary antibodies nanoparticles bind to the freely accessible Fc_γ-fragment of the user's primary antibody. The resulting hybrid forms a signaling unit that accumulates on the target protein resulting in a permanently visible signal.

Content:

Sufficient for immunostaining of five membranes, max. 100 cm² each. The primary antibody specific for the target protein is not supplied. The staining solution PIA-PINK-Rabbit is composed of anti-rabbit IgG antibody (Fc fragment, polyclonal, affinity purified, specific against the Fc_γ fragment) conjugated with gold nanoparticles (Abs_{max} 540 nm).

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Additional Storage Conditions: stable at ambient temperature for up to one week

do not freeze

Shelf Life: 6 months

Purity: > 95 % (SDS-PAGE)

Form: liquid