PDIA6 (human) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-210-889-R100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Protein disulfide-isomerase A6, Protein disulfide-

isomerase P5

Application Flow Cytometry, WB

Application NotesDetects a band of ~50kDa by Western blot.

Formulation Liquid. Neat serum containing 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 476-492

(D476IDLSDVELDDLGKDEL492) of human PDIA6 (protein disulfide-isomerase A6; protein disulfide

isomerase P5).

Recommendation Dilutions/ConditionsWestern Blot (1:1,000)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID Q15084

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

