

Lipocalin-2 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Lipocalin-2

Instruction Manual

Catalog Number	PK-AB577-3819
Description	Lipocalin 2 is from a family of proteins that are involved in the transportation of small hydrophobic molecules including steroids, retinoids, bilin and retinods. Lipocalins have been linked to many biochemical processes such as immune response, pheromone transport, biological prostaglandin synthesis, retinoid binding, and cancer cell interactions.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide corresponding to residues surrounding amino acids 40 of mouse Lipocalin 2.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-Lipocalin 2 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far. The antibody recognizes ~ 25 kDa Lipocalin 2 of human, mouse, and rat origins. Reactivity to other species has not been tested. A ~70 kDa unknown band can also be detected. Blocking peptide (Cat.# PK-AB577-3819P) is available separately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.