

Epithelial sodium channel (γ -subunit, neoepitope) monoclonal antibody (11-35-1)

The epithelial sodium channel (ENaC) of the kidney is necessary for extracellular volume homeostasis and normal arterial BP. Activity of ENaC is enhanced by proteolytic cleavage of the γ -subunit. In the γ -subunit, putative cleavage sites for prolasin, kallikrein (similar site as prolasin), elastase, and plasmin exist. Prolasin is a glycosylphosphatidylinositol anchored apical serine protease expressed in the collecting duct that activates γ ENaC at a site close to but in the C-terminal direction to the furin cleavage with putative release of a 43-residue inhibitory tract.

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

Ordering Information

[Order Online »](#)

BPD-CAM-035-01-1	1mg
BPD-CAM-035-01-04	400 μ g

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name γ ENaC

Application ELISA, IF, IHC (FS), IHC (PS), WB

Clone 11-35-1

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic peptide sequence: VGGSIIHKAC

Isotype IgG1 κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Specific for the neoepitope generated after proteolytic cleavage of human γ ENaC and recognizes the C-terminal fragment.

UniProt ID P51170

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



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)