Epithelial sodium channel (γ-subunit) monoclonal antibody (3c7)

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 and putative release of a 43-amino acid inhibitory tract from the γ -subunit ectodomain.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-CAM-003-07-1	1mg
BPD-CAM-003-07-04	400µg

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 γENaC, Amiloride-sensitive sodium channel subunit γ

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

Clone 3c7

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen The inhibitorγ peptide from the human γENaC subunit.

EAESWNSVSEGKQPRFSHRIPLC corresponding to amino acid residue 139-160 of human yENaC subunit.

Isotype IgG1ĸ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

Specificity Specific for the inhibitory tract of human γENaC subunit.

UniProt ID P51170

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

