

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](#).

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 inhibitory peptide from the human γENaC subunit. EAESWNSVSEGGKQPRFSHRIPLC corresponding to amino acid residue 139-160 of human γENaC subunit.
Isotype	IgG1κ
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](#)