CD47 recombinant monoclonal antibody (CD47/6362R)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

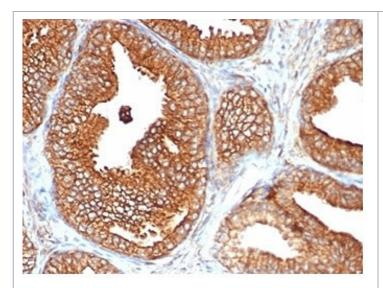
Order Online »

ENZ-ABS880-0100

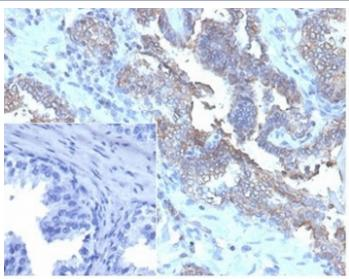
100µg

Manuals, SDS & CofA

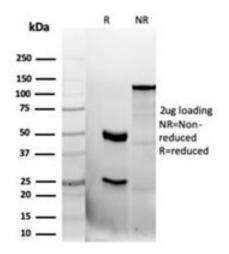
View Online »



IHC staining of FFPE human prostate carcinoma tissue with recombinant CD47 antibody (clone CD47/6362R). Negative control inset: PBS used instead of primary antibody to control for secondary Ab binding. HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human prostate carcinoma tissue with recombinant CD47 antibody (clone CD47/6362R). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



SDS-PAGE analysis of purified, BSA-free recombinant CD47 antibody (clone CD47/6362R) as confirmation of integrity and purity.

Handling & Storage

Handling Aliquot the recombinant CD47 antibody and store frozen at -20°C or colder. Avoid

freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IAP, Integrin Associated Protein

Application ELISA, IHC (PS)

Clone CD47/6362R

Formulation Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA

(US sourced), 0.05% sodium azide.

Host Rabbit

Immunogen Human CD47 (aa 18-135).

lsotype lgGk

Purity Detail Protein A/G affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (FFPE) (1-2µg/ml)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID Q08722

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

•

Last modified: July 12, 2024



info-

eu@enzolifesciences.com