CD44v4 Recombinant monoclonal antibody (CD44v4/1700R)

The CD44 molecule belongs to a family of cellular adhesion molecules found on a wide range of normal and malignant cells in epithelial, mesothelial and hemopoiesis tissues. It is a single gene with 20 exons, of which 10 are normally expressed to encode the basic CD44 (H-CAM) molecule. The additional 10 exons (v1 to v10) are only expressed by alternative splicing of the nuclear RNA. The expression of specific cell adhesion molecule CD44 splice variants has been reported to be associated with metastasis in certain human malignancies.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS349-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, IF, IHC (PS)

Clone CD44v4/1700R

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced)

and 0.05% sodium azide.

Host Rabbit

Immunogen An amino acid sequence from the variant 4 domain of

CD44.

Isotype IgGk

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (0.5-1 μg/106 cells in 0.1

mL)Immunofluorescence (0.5-1

 $\mu g/mL) lmmunohistochemistry (formalin-fixed paraffin sections) (0.5-1 <math display="inline">\mu g/mL$ for 30 min at RT)Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

SpecificityThis antibody recognizes an epitope encoded by exon v4

on the variant portion of human CD44.

UniProt ID P16070

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com