CD40 (human) monoclonal antibody (G28-5)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

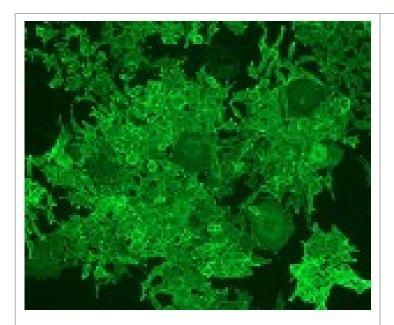
Order Online »

ENZ-ABS130-0100

100µg

Manuals, SDS & CofA

View Online »



Immunohistochemistry analysis of CD40-transfected L cells stained with CD40 (human) mAb, clone G28-5.

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tumor necrosis factor receptor superfamily member 5,

TNFRSF 5, B-cell surface antigen CD40, Bp50, CD40L

receptor, CDw40

Application Flow Cytometry, FUNC, IHC (PS)

Application Notes Functional Application: cell activation.

Clone G28-5

Formulation Liquid. In TRIS-NaCl, pH 8.

Host Mouse

Immunogen Human erythrocyte-rosette-negative tonsillar lymphocyte.

Isotype IgG1k

Purity Detail Purified using QMA Hyper D ion exchange

chromatography

Species Reactivity Human

UniProt ID P25942

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com