



PRODUCT DATA SHEET

CD30 Recombinant monoclonal antibody (Ki-1/1505R)

ENZ-ABS354

Product Number/Sizes

ENZ-ABS354-0100 100 µg

Product Details

ALTERNATIVE NAME: Ki-1, TNFRSF 8
CLONE: Ki-1/1505R
HOST: Rabbit
ISOTYPE: IgGκ
IMMUNOGEN: Partial human CD30.
UNIPROT ID: P28908
SPECIES REACTIVITY: Human
APPLICATIONS: Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS: Flow Cytometry (0.5-1 µg/10⁶ cells in 0.1 mL)
Immunofluorescence (1-2 µg/mL)
Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/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.
PURITY DETAIL: Protein A affinity purified.
FORMULATION: Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
SHIPPING: Blue Ice
LONG TERM STORAGE: +4°C
SCIENTIFIC BACKGROUND: CD30 is the receptor for TNFSF8/CD30L. May play a role in the regulation of cellular growth and transformation of activated lymphoblasts. Regulates gene expression through activation of NF-kappa-B. [UniProt]
REGULATORY STATUS: RUO - Research Use Only

Revised 19-Dec-23

GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.
10 Executive Blvd
Farmingdale, NY 11735
USA
T 1-800-942-0430
T 1-631-694-7070
F 1-610-941-9252
E info-usa@enzolifesciences.com
www.enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lausen
Switzerland
T +41/061 926 89 89
F +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com