



# PRODUCT DATA SHEET

## CD30 Recombinant monoclonal antibody (Ki-1/1747R)

ENZ-ABS355

### Product Number/Sizes

ENZ-ABS355-0100 100 µg

### Product Details

ALTERNATIVE NAME: Ki-1, TNFRSF 8  
CLONE: Ki-1/1747R  
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/10<sup>6</sup> 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