

CD30 Recombinant monoclonal antibody (rKi-1/779)

ENZ-ABS327

Product Number/Sizes

ENZ-ABS327-0100 100 µg

Product Details

ALTERNATIVE NAME: Ki-1, TNFRSF 8
CLONE: rKi-1/779
HOST: Mouse
ISOTYPE: IgG1κ
IMMUNOGEN: Recombinant full length 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.
APPLICATION NOTES: This mAb distinguishes large cell lymphomas derived from activated lymphoid cells from histiocytic malignancies and lymphomas derived from resting and precursor lymphoid cells or from anaplastic carcinomas.
PURITY DETAIL: Protein G 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 synthesized as a 90kDa precursor, which is processed in the Golgi complex into a membrane-bound phosphorylated mature 105/120kDa glycoprotein. In Hodgkin's disease, CD30/Ki-1 antigen is expressed by mononuclear-Hodgkin and multinucleated Reed-Sternberg cells. It is also expressed by the tumor cells of a majority of anaplastic large cell lymphomas as well as by a varying proportion of activated T and B cells. About one third of the CD30/Ki-1 positive lymphomas lack the leukocyte common antigen (CD45).
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