



PRODUCT DATA SHEET

CA19-9 Recombinant monoclonal antibody (CA19.9/1390R)

ENZ-ABS356

Product Number/Sizes

ENZ-ABS356-0100 100 µg

Product Details

CLONE:	CA19.9/1390R
HOST:	Rabbit
ISOTYPE:	IgGκ
IMMUNOGEN:	Purified human CA19-9.
SPECIES REACTIVITY:	Human
APPLICATIONS:	Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS:	Flow Cytometry (2-4 µg/10 ⁶ cells in 0.1 mL) Immunofluorescence (5-10 µg/mL) Immunohistochemistry (formalin-fixed paraffin sections) (5-10 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:	CA19-9, a carbohydrate epitope expressed on a high MW (>400kDa) mucin glycoprotein, is a sialyl Lewis a structure which is synthesized from type 1 blood group precursor chains and is present in individuals expressing the Lewis a and/or Lewis b blood group antigens.
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