

CD303 monoclonal antibody (124B3.13) (purified)

NOT SOLD IN GERMANY.

Ordering Information

[Order Online »](#)

DEN-DDX0043P-100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name BDCA-2, Blood dendritic cell antigen 2

Application IHC (PS)

Application Notes Immunohistochemistry (formol paraffine)

Clone 124B3.13

Host Mouse

Isotype IgG1

Species Reactivity Human, Mouse, Rat

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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