

DC-LAMP/CD208 monoclonal antibody (1010E1.01) (purified)

Ordering Information

[Order Online »](#)

DEN-DDX0191P-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

Application Flow Cytometry, IHC (PS)

Application Notes Immunohistochemistry (formol paraffine, bouin paraffine), Flow Cytometry (intracyto in human)

Clone 1010E1.01

Host Rat

Isotype IgG2a

Species Reactivity Cat, Dog, Human, Mouse, Sheep

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