

DLC1 monoclonal antibody (10D6)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-3400	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Dynein light chain 1, PIN, Protein inhibitor of nNOS, LC8
Application	Flow Cytometry, WB
Clone	10D6
Crossreactivity	Does not cross-react with DLC2.
Formulation	Liquid. In PBS containing 10mM sodium azide.
Host	Rat
Immunogen	Recombinant mouse DLC1 (dynein light chain 1; LC8; PIN).
Isotype	IgM
Purity Detail	Affinity purified.
Species Reactivity	Human, Mouse
UniProt ID	P63168
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



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

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

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

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

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

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

Last modified: May 29, 2024