

DLC1 monoclonal antibody (10D6)

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-804-340-C100 | 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](#)

.



Last modified: May 29, 2024

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
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420

info@enzolifesciences.com
be@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
uk@enzolifesciences.com