

Msx2 monoclonal antibody (2E12)

Msx2, also known as HOX-8, is a homeobox containing transcription factor, which is a member of the MSH family of homeobox proteins. Homeobox (HOX) genes encode DNA binding proteins that recognize specific sequences and modulate transcriptional activity. Msx2 is believed to play a role in regulating bone development, particularly in sutures of cranial bone.

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

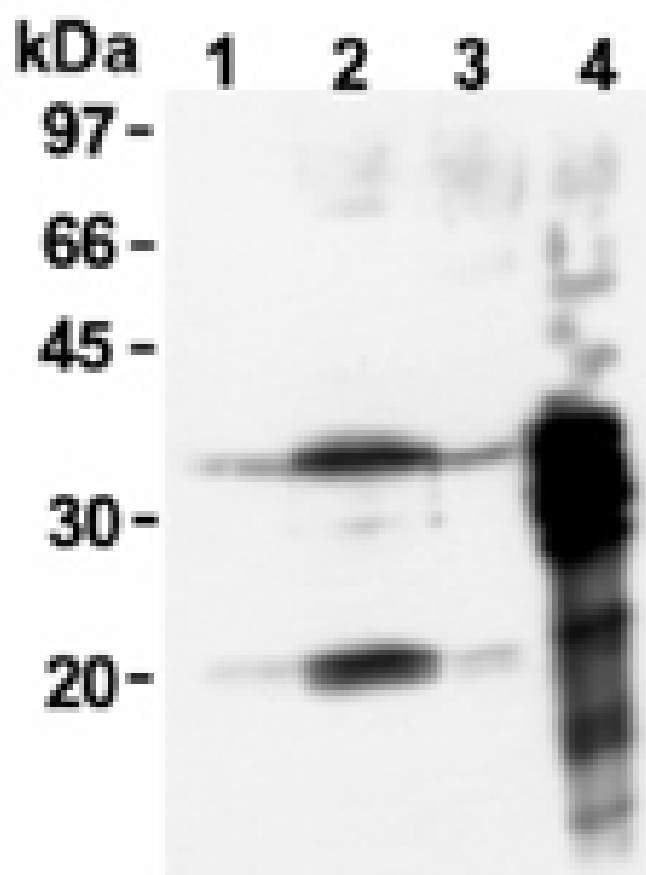
Ordering Information

[Order Online »](#)

ADI-AAM-010-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HeLa (1), ZR-75-1 (2), KB (3) cell lysates, and recombinant Msx2 protein (4), probed with Msx2 mAb.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Homeobox protein MSX-2, Hox-8
Application	IHC, WB
Application Notes	Detects a band of ~30kDa by Western blot.
Clone	2E12
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
Host	Mouse
Immunogen	Recombinant human Msx2 (N-terminal fragment).
Isotype	IgG2a
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (10µg/ml)Western Blot (1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P35548
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
usa@enzolifesciences.com

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