

Polycomb 2

polyclonal antibody

Polycomb 2 is a member of *Drosophila* Polycomb group gene family. The polycomb group (PcG) genes are essential for maintenance of proper expression patterns of developmental master regulators; changes in expression of PcG proteins have been associated with cancer. Polycomb 2 is a part of the cellular memory system responsible for the inheritance of gene activity by progeny cells. It participates in maintaining the transcriptionally repressive state of genes. Polycomb 2 is part of a complex that acts via chromatin remodeling and modification of histones; it mediates monoubiquitination of histone H2A 'Lys-119', rendering chromatin heritably changed in its expressibility. Polycomb 2 is an E3 SUMO-protein ligase which facilitates SUMO1 conjugation by UBE2I.

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

Ordering Information

[Order Online »](#)

ENZ-ABS291-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Aliquots are stable for at least 12 months.
Handling	Avoid freeze/thaw cycles, aliquot and store at -20°C. For maximum recovery of product, centrifuge the original vial after thawing and prior to removing the cap.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	E3 SUMO-protein Ligase CBX4, Chromobox Protein Homolog 4, Pc2
Application	ELISA, IHC, WB
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of N-terminal human Polycomb 2, conjugated to KLH.
Isotype	IgG
Purity Detail	Affinity purified
Recommendation Dilutions/Conditions	ELISA (1:1,000)Immunohistochemistry (1:50-1:100)Western Blot (1:100-1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	O00257

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-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