

Phosphoinositide 3-kinase p110 δ (human) polyclonal antibody

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

Ordering Information

[Order Online »](#)

ALX-210-747-C200	200 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	PI(3)K p110δ
Application	ICC, IHC (PS), WB
Application Notes	Detects a band of ~110kDa by Western blot.
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 1026-1044 (S1026WKTKVNWLAHNVS KDNRQ1044) of the C-terminal domain of human PI(3)K p110δ (phosphoinositide 3-kinase p110δ).
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Western blot on human T cell line Jurkat.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:300)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	O00329

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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