

# Phosphoinositide 3-kinase p110 $\delta$ (human) polyclonal antibody

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

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Handling</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	PI(3)K p110δ
<b>Application</b>	ICC, IHC (PS), WB
<b>Application Notes</b>	Detects a band of ~110kDa by Western blot.
<b>Formulation</b>	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
<b>Host</b>	Rabbit
<b>Immunogen</b>	Synthetic peptide corresponding to aa 1026-1044 (S1026WKTKVNWLAHNVSKDNRQ1044) of the C-terminal domain of human PI(3)K p110δ (phosphoinositide 3-kinase p110δ).
<b>Purity Detail</b>	Epitope-affinity purified IgG.
<b>Quality Control</b>	Western blot on human T cell line Jurkat.
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	O00329
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)