

PIDD monoclonal antibody (Anto-1)

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

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-837-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

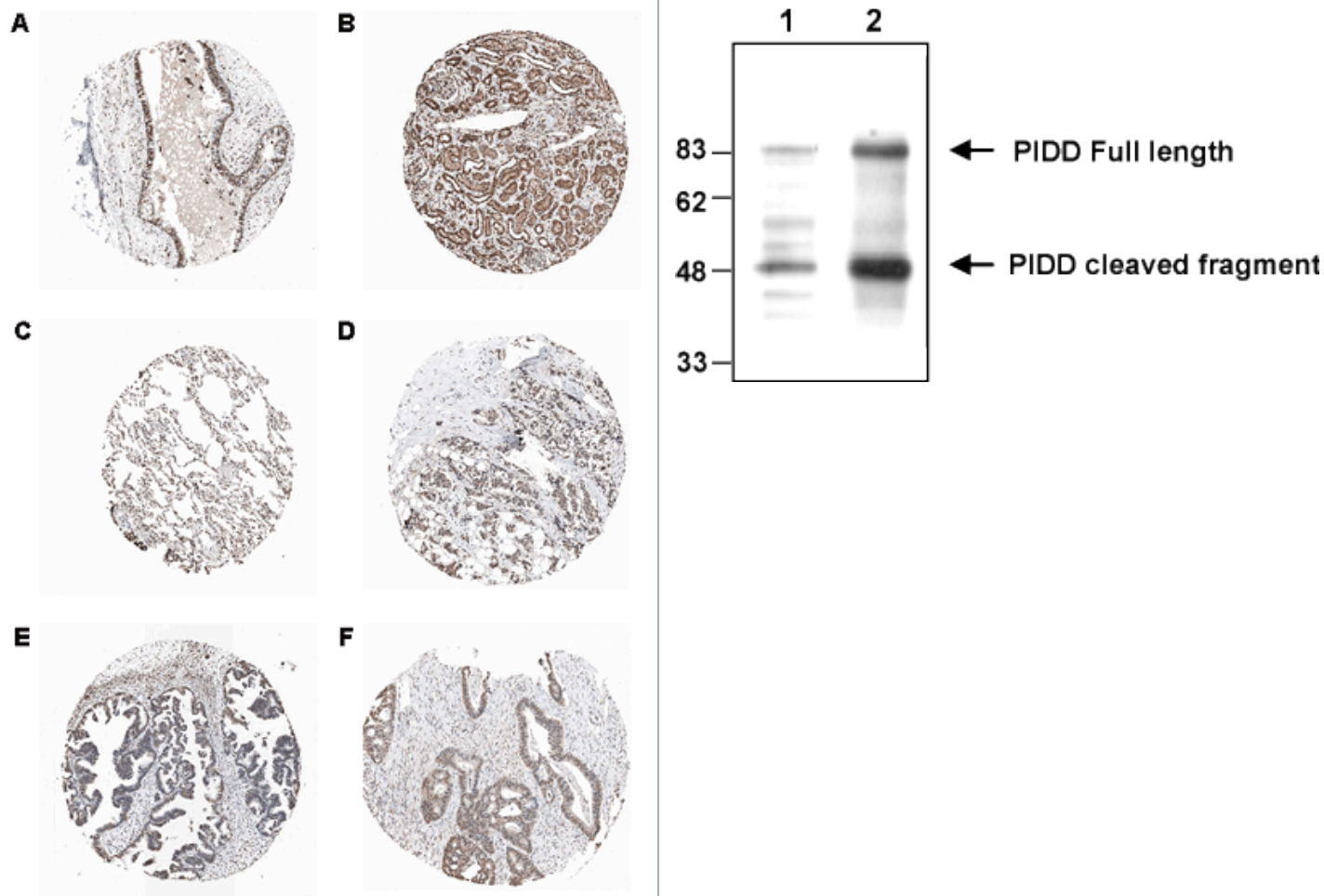


Figure 2: Immunohistochemical staining of endogenous human PIDD in different human tissues (paraffin sections) using MAb to PIDD (Anto-1) (Prod. No. [ALX-804-837](#)). **Method:** Different human normal (A: Bronchus; B: Kidney; C: Lung) or cancer (D: Breast; E: Ovarian; F: Colo-rectal) tissues were stained with MAb to PIDD (Anto-1) by standard immunohistochemistry. *Pictures courtesy of Human Protein Atlas (*www.proteinatlas.org*).* “””””

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IHC (PS), WB
Application Notes	Detects bands of ~90kDa (full length protein) and ~50kDa (cleaved protein) by Western blot.
Clone	Anto-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human PIDD death domain (aa 776-910).
Isotype	IgG1
Recommendation Dilutions/Conditions	Western Blot (1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Human, Mouse
UniProt ID	Q9HB75



ENZO LIFE SCIENCES,
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 1 402 44 655
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

Worry Free Guarantee This antibody is covered by our **Worry Free Guarantee**