

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 »](#)

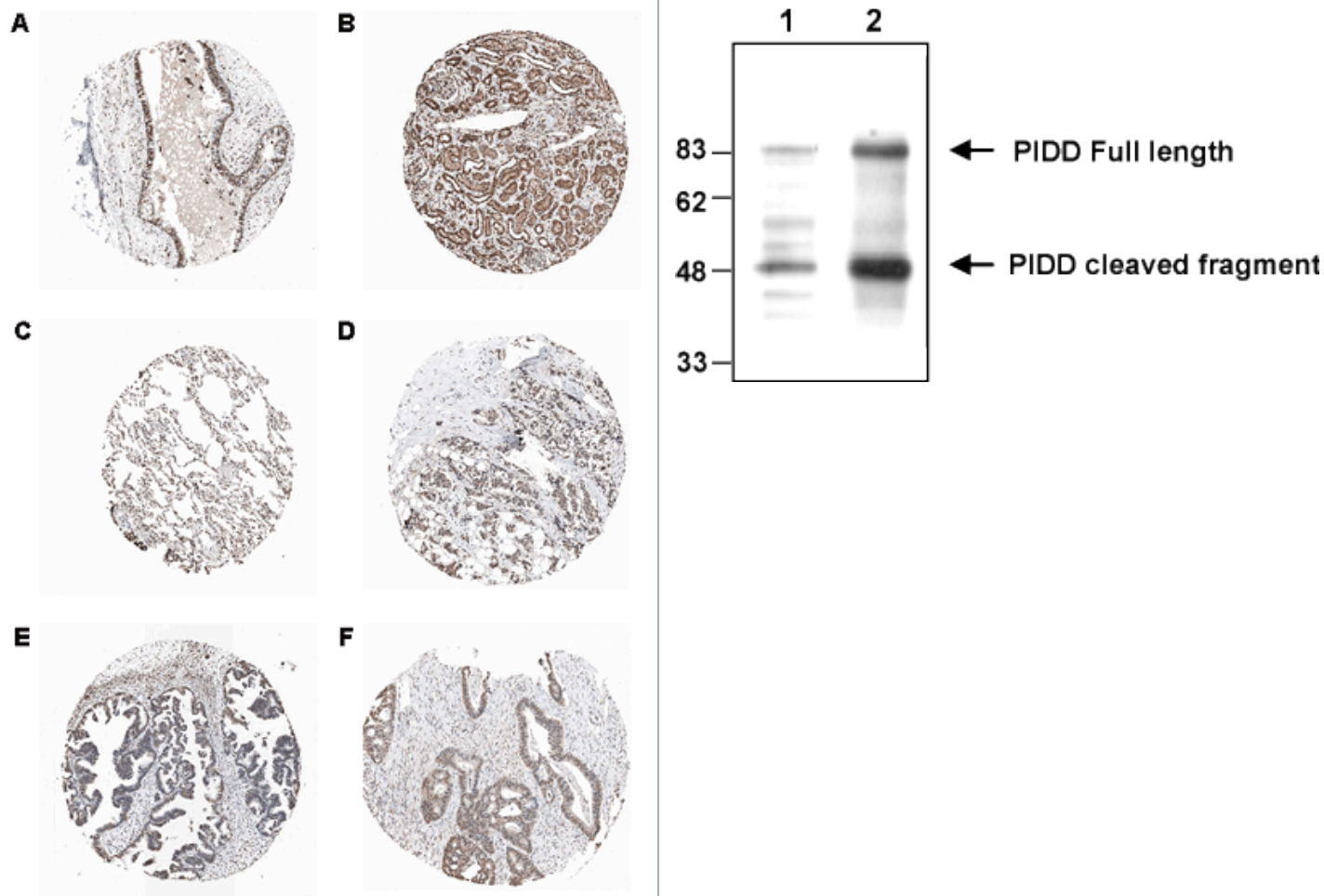


Figure 2: Immunohistochemical staining of endogenous human PIDD in different human tissues (paraffin sections) using MAb to PIDD (Anto-1) (Prod. No. [ALX-804-837](#)). **Method:** Different human normal (A: Bronchus; B: Kidney; C: Lung) or cancer (D: Breast; E: Ovarian; F: Colo-rectal) tissues were stained 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
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)