

Placental protein 14 (human) monoclonal antibody (16)

Human placental protein 14 (PP14; also glycodelin and progesterone-associated endometrial protein) is a protein of the lipocalin structural superfamily. PP14 is the most abundant product of the secretory endometrium, and has been proposed as a biochemical marker of endometrial function in women undergoing in vitro fertilization treatment.

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

Ordering Information

[Order Online »](#)

BPD-BTE-001-16-1	1mg
BPD-BTE-001-16-04	400µg

Manuals, SDS & CofA

[View Online »](#)

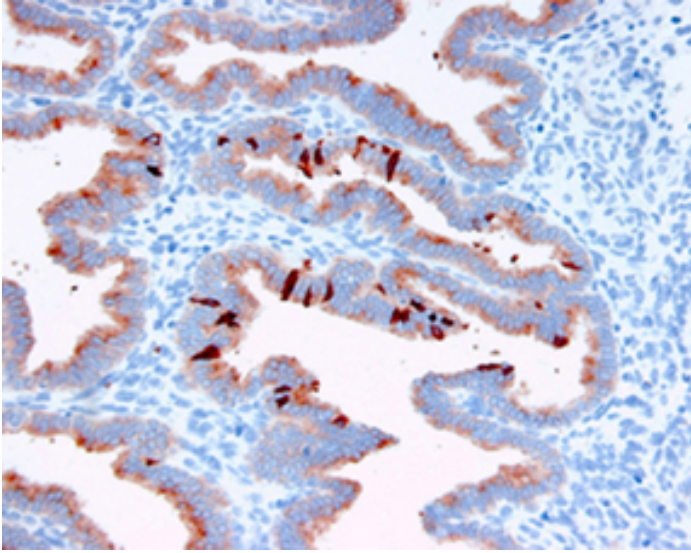


Figure: Immunohistochemical analysis of formalin-fixed, paraffin-embedded late secretory phase endometrium tissue using Placental protein 14 (human), mAb. Specific staining (red-brown) is restricted to the glandular cells of the endometrium.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Glycodelin A, GD, Pregnancy-associated endometrial alpha-2 globulin, PEG, PAEG, PP14, Progesterone-associated endometrial protein, Progestagen-associated endometrial protein
Application	ELISA, IHC
Clone	16
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	5047
Host	Mouse
Immunogen	Purified human placental protein 14 (glycodelin A) from second trimester amniotic fluid.
Isotype	IgG1κ
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
UniProt ID	P09466

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)