

# Placental protein 14 (human) monoclonal antibody (13)

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-001-13-1	1mg
BPD-BTE-001-13-04	400µg

## Manuals, SDS & CofA

[View Online »](#)

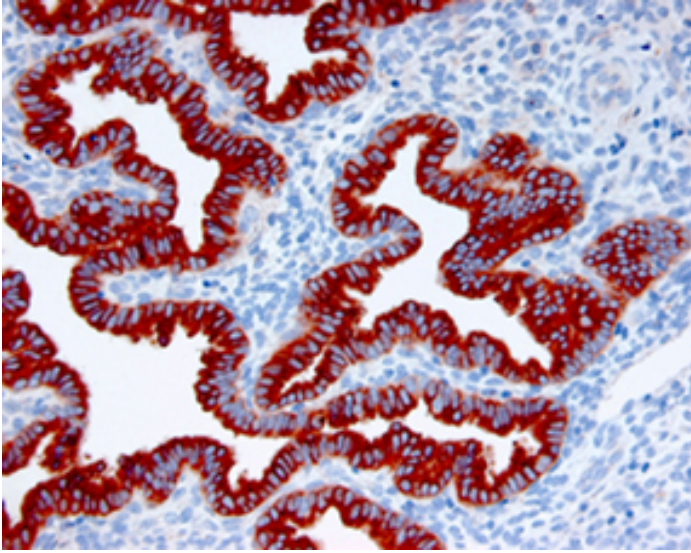


Figure: Immunohistochemical analysis of formalin-fixed paraffin-embedded endometrial tissue from the late secretory phase using Placental protein 14 (human), mAb. Staining is restricted to the glandular cells.

## 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	13
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.
Species Reactivity	Human
UniProt ID	P09466

## Worry-free Guarantee

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



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 940 946 9120  
[usa@enzolifesciences.com](mailto: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)