

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](#).

Manuals, SDS & CofA

[View Online »](#)

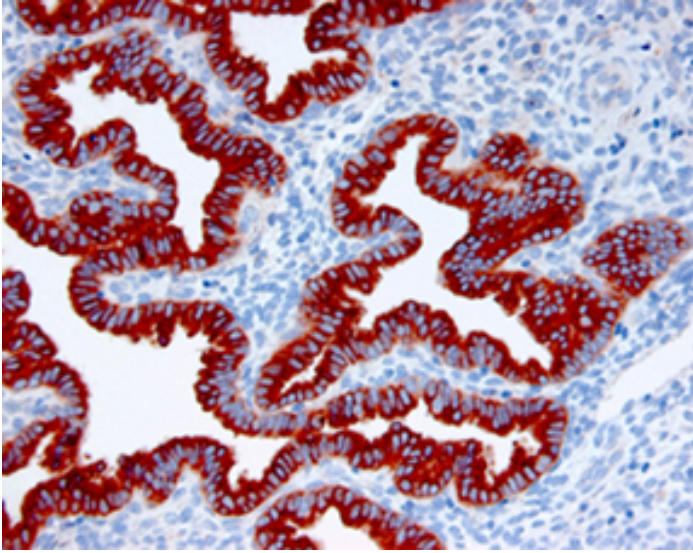


Figure: Immunohistochemical analysis of formalin-fixed paraffin-embedded endometrial tissue from the late secretory phase using Placental protein 14 (human), mAb. Staining is restricted to the glandular cells.

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	IgG1k
Purity Detail	Protein A/G-purified.
Species Reactivity	Human
UniProt ID	P09466
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee