



Product datasheet

Human FOLR3 ELISA Kit ab267613

1 References 1 Image

Overview

Product name	Human FOLR3 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Serum, Plasma, Cell culture media				
Assay type	Sandwich (quantitative)				
Sensitivity	32 pg/ml				
Range	32.77 pg/ml - 8000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		103		96% - 110%
	Plasma		96		86% - 108%
	Cell culture media		91		89% - 95%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human FOLR3 ELISA Kit is designed for the quantitative determination of FOLR3 in cell culture supernatants, plasma and serum samples.				

This assay employs an antibody specific for human FOLR3 coated on a 96-well plate. Standards and samples are pipetted into the wells and FOLR3 present in a sample is bound to the wells by

the immobilized antibody. The wells are washed and biotinylated anti-human FOLR3 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of FOLR3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Anti-Human FOLR3 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human FOLR3 Detection Antibody	2 vials
Human FOLR3 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

Function Binds to folate and reduced folic acid derivatives and mediates delivery of 5-methyltetrahydrofolate to the interior of cells. Isoform Short does not bind folate.

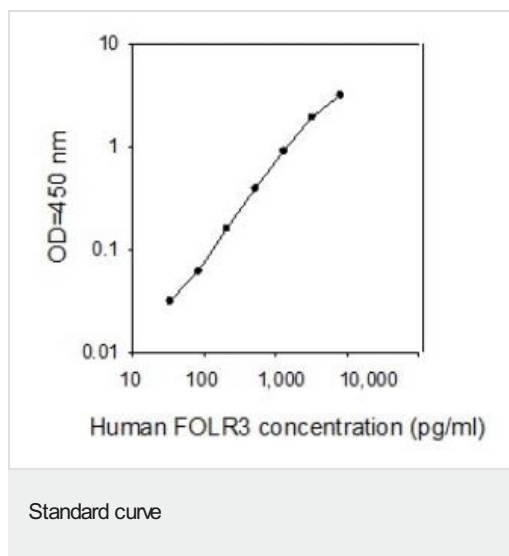
Tissue specificity Spleen, thymus, bone marrow, ovarian carcinoma, and uterine carcinoma.

Sequence similarities Belongs to the folate receptor family.

Post-translational modifications Eight disulfide bonds are present.

Cellular localization Secreted.

Images



Human FOLR3 ELISA kit (ab267613) Standard curve.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors