

Rat GPNMB ELISA Kit ab270219

リコンビナント SimpleStep ELISA

画像数 5

製品の概要

製品名 Rat GPNMB ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Serum	7			4.3%

サンプルの種類 Cell culture supernatant, Serum, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 136 pg/ml

検出範囲 0.31 ng/ml - 20 ng/ml

添加回収試験 特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	104	94% - 122%
Serum	98	95% - 101%
Cell culture media	112	109% - 114%
Hep Plasma	105	94% - 121%
EDTA Plasma	102	97% - 112%
Cit plasma	120	112% - 133%

全工程の試験時間 1h 30m

ステップ One step assay

種交差性 交差種: Rat

製品の概要 Rat GPNMB ELISA Kit (ab270219) is a single-wash 90 min sandwich ELISA designed for the

quantitative measurement of GPNMB protein in cell culture media, cell culture supernatant, edta plasma, serum, cit plasma, and hep plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Rat GPNMB with 136 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

特記事項

GPNMB (Glycoprotein non-metastatic gene B), also known as Osteoactivin, is a type I transmembrane glycoprotein most often associated with endosomal/lysosomal compartment and over expressed in many types of cancer. The GPNMB is involved in many physiological processes, including transport of late melanosomes to keratinocytes, regulation of osteoblast and osteoclast, stimulation of dendritic cell maturation, and stimulation of adhesion of dendritic cells to endothelial cells.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

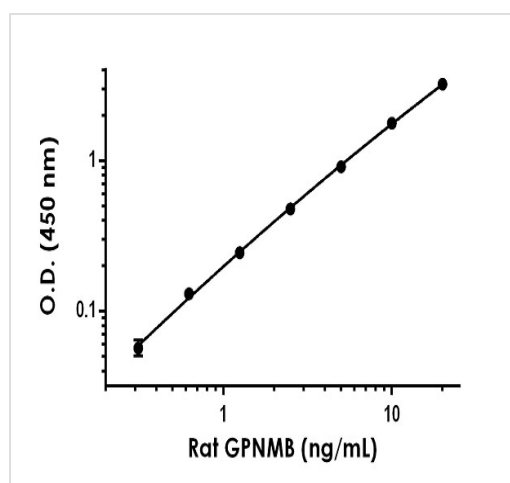
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Rat GPNMB Capture Antibody	1 x 600µl
10X Rat GPNMB Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Plate Seals	1 unit
Rat GPNMB Lyophilized Recombinant Protein	2 vials
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml

内容	1 x 96 tests
TMB Development Solution	1 x 12ml

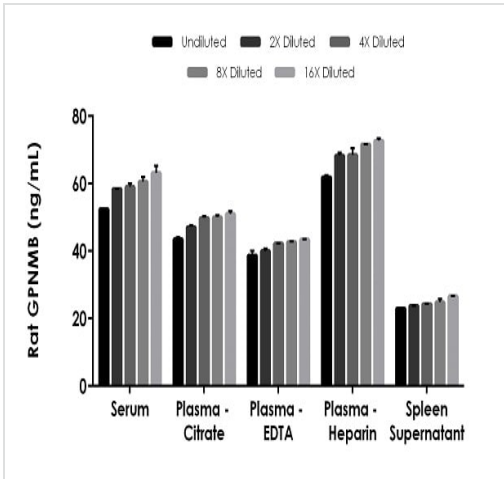
機能	Could be a melanogenic enzyme.
組織特異性	Not restricted to the melanocytic lineage.
配列類似性	Belongs to the PMEL/NMB family. Contains 1 PKD domain.
発生段階	Expression in poorly metastatic melanoma cell lines; no expression in highly metastatic melanoma cell lines.
細胞内局在	Membrane. Melanosome. Identified by mass spectrometry in melanosome fractions from stage I to stage IV.

画像



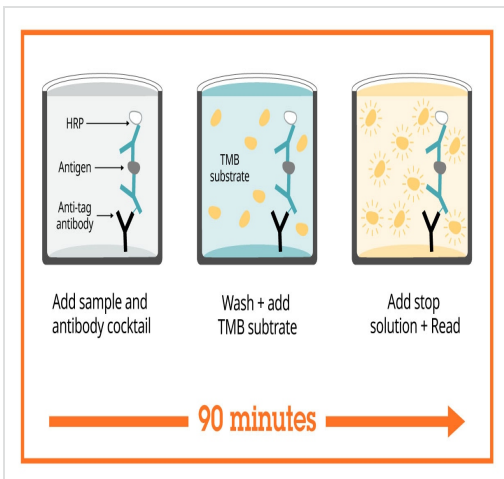
The GPNMB standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of rat GPNMB standard curve in Sample Diluent NS.



Interpolated concentrations of native GPNMB in rat serum, plasma and cell culture supernatant samples.

The concentrations of GPNMB were measured in duplicates, interpolated from the GPNMB standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 20%, plasma (citrate) 20%, plasma (EDTA) 40%, plasma (heparin) 20% and spleen supernatant 80%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean GPNMB concentration was determined to be 59 ng/mL in neat serum, 48 ng/mL in neat plasma (citrate), 41 ng/mL in neat plasma (EDTA), 69 ng/mL in neat plasma (heparin) and 24 ng/mL in neat spleen supernatant.



Sandwich ELISA - Rat GPNMB ELISA Kit (ab270219)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Powered by recombinant antibodies

- Research with confidence**
Consistent and reproducible results
- Long-term and scalable supply**
Recombinant technology
- Success from the first experiment**
Confirmed specificity
- Ethical standards compliant**
Animal-free production

Sandwich ELISA - Rat GPNMB ELISA Kit (ab270219)

To learn more about the advantages of recombinant antibodies see [here](#).

Get more done with SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Rat GPNMB ELISA Kit
(ab270219)

To learn more about the advantages of SimpleStep ELISA® kits see [here](#).

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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