

Serotonin ELISA Kit ab133053

★★★★☆ 4 Abreviews 28 References 画像数 3

製品の概要

| | |
|------|---------------------|
| 製品名 | Serotonin 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% |
|-------------|----|-----|----|------|
| 346.1 ng/ml | 20 | | | 4.2% |
| 53.8 ng/ml | 20 | | | 5.8% |
| 13.5 ng/ml | 20 | | | 11% |

Inter-Assay (日差再現性)

| サンプル | N | 平均値 | SD | CV% |
|-------------|---|-----|----|-------|
| 358.0 ng/ml | | | | 16.2% |
| 53.9 ng/ml | | | | 18.4% |
| 13.2 ng/ml | | | | 12.7% |

| | |
|---------|---------------------------------|
| サンプルの種類 | Urine, Serum, Plasma, Platelets |
| アッセイタイプ | Competitive |
| 検出感度 | 0.293 ng/ml |
| 検出範囲 | 0.49 ng/ml - 500 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 |
|-----------|------|-------------------|
| Urine | 101 | 5ng/ml - 100ng/ml |
| Serum | 91 | 5ng/ml - 100ng/ml |
| Platelets | 107 | 5ng/ml - 100ng/ml |

| サンプルの種類 | 平均 % | 測定範囲 |
|------------|------|-------------------|
| Cit plasma | 95 | 5ng/ml - 100ng/ml |

全工程の試験時間

3h 00m

ステップ

Multiple steps standard assay

種交差性

交差種: Species independent

製品の概要

Serotonin ELISA kit is a competitive ELISA kit designed for the accurate quantitative measurement of Serotonin in Human Serum, Platelets, Plasma and Urine.

A goat anti-rabbit IgG antibody has been precoated onto 96-well plates. Standards or test samples are added to the wells, along with an alkaline phosphatase (AP) conjugated- Serotonin antigen and a polyclonal rabbit antibody specific to Serotonin. After incubation the excess reagents are washed away. pNpp substrate is added and after a short incubation the enzyme reaction is stopped and the yellow color generated is read at 405 nm. The intensity of the yellow coloration is inversely proportional to the amount of Serotonin captured in the plate.

特記事項

Serotonin (5-hydroxytryptamine, 5-HT) is a monoamine found in the central nervous system, gastrointestinal tract, and blood with broad physiological functions as a neurotransmitter, in gastric motility, hemostasis, and cardiovascular integrity. Defects in serotonin signaling have been linked to a large number of complex behavioral disorders in humans, including anxiety and depression, autism, and eating disorders, and to neurodegenerative disorders such as Parkinson's. Platelets serve as the major reservoir of serotonin in the bloodstream. When activated, platelets release serotonin into the bloodstream where it acts as a powerful vasoconstrictor. Treatment with Selective Serotonin Uptake Inhibitors (SSRIs) dramatically reduces platelet serotonin concentrations, and altered levels of serotonin in the circulatory system are implicated in such diverse conditions such as asthma, liver cirrhosis and carcinoid tumors.

Cross reactivities:

| Compound | Cross Reactivity |
|----------------------------|------------------|
| N-Acetyl Serotonin | 17% |
| 5-Hydroxy-L-tryptophan | 0.4% |
| Tryptamine | 0.1% |
| 5-Hydroxyindoleacetic acid | 0.03% |
| Melatonin | 0.01% |
| Tyramine | <0.004% |
| Tryptophan | <0.004% |

試験プラットフォーム

Microplate

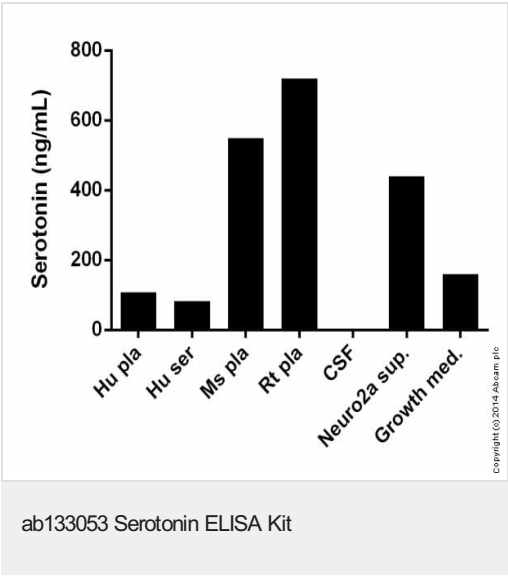
製品の特長

保存方法

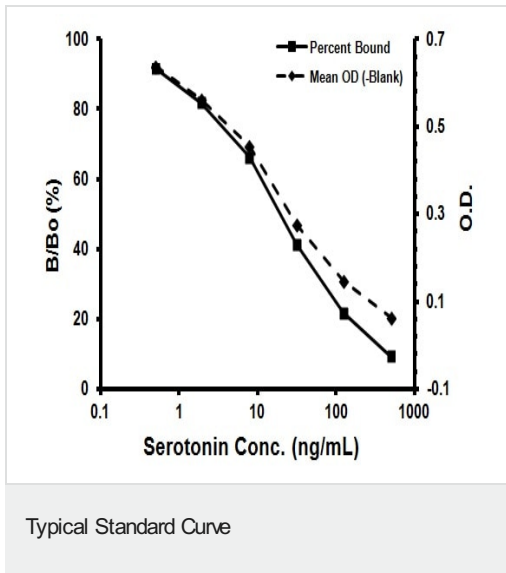
Store at -20°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer Concentrate | 1 x 27ml |
| Assay Buffer | 1 x 27ml |
| Goat anti-rabbit IgG Microplate (12 x 8 wells) | 1 unit |
| Plate Sealer | 2 units |
| pNpp Substrate | 1 x 20ml |
| Serotonin Alkaline Phosphatase Conjugate | 2 x 150ng |
| Serotonin Antibody | 2 vials |
| Serotonin Standard | 2 x 250ng |
| Stop Solution | 1 x 5ml |

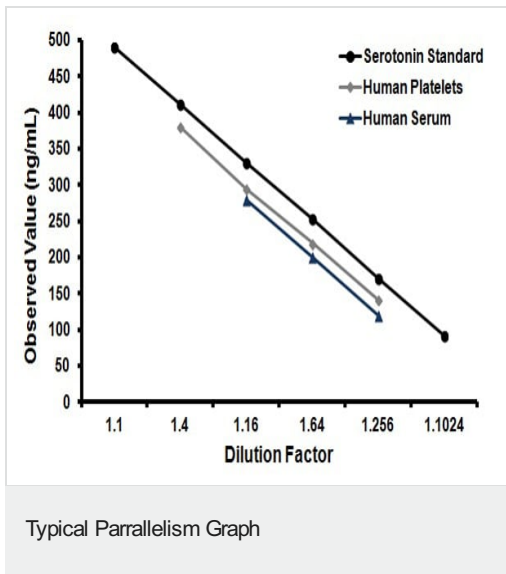
画像



Serotonin measured (averages of undil., 1:3, 1:9 & 1:27) in biological fluids (Neuro2a an Growth med. averaged from 1:2 and 1:4) showing quantity (ng) per mL of tested sample.



Representative Standard Curve using ab133053.



Dose-response curves from Human platelets and Human serum diluted into assay buffer were compared to the Serotonin standard curve. The parallel response indicates the standard effectively mimics the native protein.

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