

Human Sonic Hedgehog ELISA Kit ab100639

★★★★★ 1 Abreviews 23 References 画像数 2

製品の概要

| | |
|---------|---|
| 製品名 | Human Sonic Hedgehog ELISA Kit |
| 検出方法 | Colorimetric |
| サンプルの種類 | Cell culture supernatant, Serum, Plasma |
| アッセイタイプ | Sandwich (quantitative) |
| 検出感度 | < 20 pg/ml |
| 検出範囲 | 13.72 pg/ml - 10000 pg/ml |
| 添加回収試験 | 94 % |

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 | 91.72 | 82% - 101% |
| Serum | 95.34 | 84% - 103% |
| Plasma | 76.43 | 68% - 89% |

| | |
|------------|---|
| ステップ | Multiple steps standard assay |
| 種交差性 | 交差種: Human |
| 製品の概要 | <p>Abcam's Sonic Hedgehog Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human Sonic Hedgehog in serum, plasma, cell culture supernatants.</p> <p>This assay employs an antibody specific for Human Sonic Hedgehog coated on a 96-well plate. Standards and samples are pipetted into the wells and Sonic Hedgehog present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Sonic Hedgehog 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 Sonic Hedgehog bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p> |
| 特記事項 | Optimization may be required with urine samples. |
| 試験プラットフォーム | Microplate |

製品の特性

保存方法

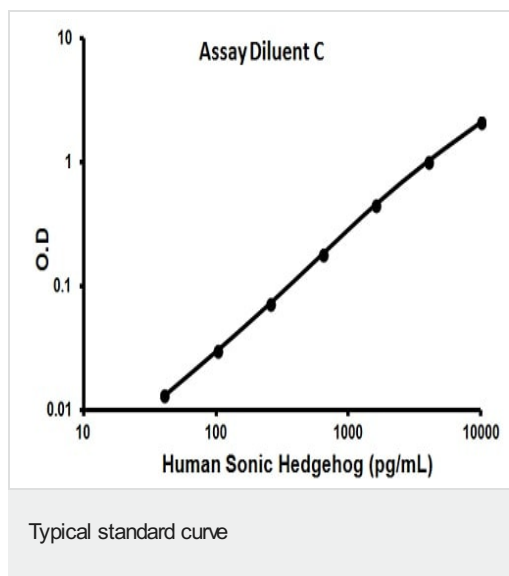
Store at -20°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|---|--------------|
| 200X HRP-Streptavidin Concentrate | 1 x 200µl |
| 20X Wash Buffer | 1 x 25ml |
| 5X Assay Diluent B | 1 x 15ml |
| Assay Diluent C | 1 x 30ml |
| Biotinylated anti-Human Sonic Hedgehog | 2 units |
| Recombinant Human Sonic Hedgehog Standard (lyophilized) | 2 vials |
| Sonic Hedgehog Microplate (12 x 8 wells) | 1 unit |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

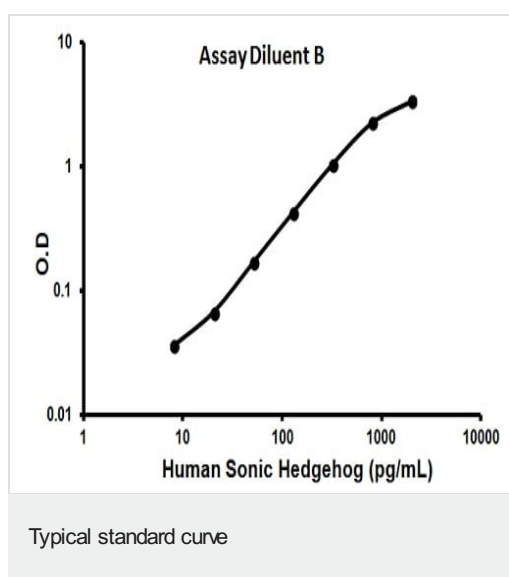
関連性

Sonic Hedgehog, which is expressed only during embryogenesis, is instrumental in patterning the early embryo. It has been implicated as the key inductive signal in patterning of the ventral neural tube, the anterior-posterior limb axis, and the ventral somites. Of three human proteins showing sequence and functional similarity to the Sonic Hedgehog protein of *Drosophila*, this protein is the most similar. The protein is made as a precursor that is autocatalytically cleaved; the N-terminal portion is soluble and contains the signalling activity while the C-terminal portion is involved in precursor processing. More importantly, the C-terminal product covalently attaches a cholesterol moiety to the N-terminal product, restricting the N-terminal product to the cell surface and preventing it from freely diffusing throughout the developing embryo. Defects in this protein or in its signalling pathway are a cause of holoprosencephaly (HPE), a disorder in which the developing forebrain fails to correctly separate into right and left hemispheres. HPE is manifested by facial deformities. In addition, it is thought that mutations in this gene or in its signalling pathway may be responsible for VACTERL syndrome, which is characterized by vertebral defects, anal atresia, tracheoesophageal fistula with esophageal atresia, radial and renal dysplasia, cardiac anomalies, and limb abnormalities.

画像



Representative standard curve using ab100639



Representative standard curve using ab100639

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