

Human LOXL2 ELISA Kit ab213808

1 References 画像数 1

製品の概要

製品名	Human LOXL2 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%
	1	16	252pg/ml	11.84	4.7%
	2	16	1465pg/ml	114.27	7.8%
	3	16	3854pg/ml	254.36	6.6%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	1	24	263pg/ml	13.93	5.3%
	2	24	1490pg/ml	123.67	8.3%
	3	24	4228pg/ml	317.1	7.5%
サンプルの種類	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 10 pg/ml				
検出範囲	156 pg/ml - 10000 pg/ml				
全工程の試験時間	3h 30m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	The Human LOXL2 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213808) is designed				

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

polyclonal antibody from goat specific for LOXL2 has been pre-coated onto 96-well plates. Standards (NSO, Q26-Q774) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for LOXL2 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human LOXL2 amount of sample captured in plate.

特記事項

Lysyl oxidase homolog 2 is an enzyme that in humans is encoded by the LOXL2 gene. This gene encodes a member of the lysyl oxidase gene family. The prototypic member of the family is essential to the biogenesis of connective tissue; encoding an extracellular copper-dependent amine oxidase that catalyzes the first step in the formation of crosslinks in collagens and elastin. A highly conserved amino acid sequence at the C-terminus end appears to be sufficient for amine oxidase activity, suggesting that each family member may retain this function. The N-terminus is poorly conserved and may impart additional roles in developmental regulation, senescence, tumor suppression, cell growth control and chemotaxis to each member of the family. LOXL2 can also crosslink collagen type IV and hence influence the sprouting of new blood vessels.

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 -20°C. Please refer to protocols.

内容	ラベル	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human LOXL2 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human LOXL2 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human LOXL2 standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

組織特異性

Expressed in many tissues. Highest expression in reproductive tissues, placenta, uterus and prostate.

配列類似性

Belongs to the lysyl oxidase family.

Contains 4 SRCR domains.

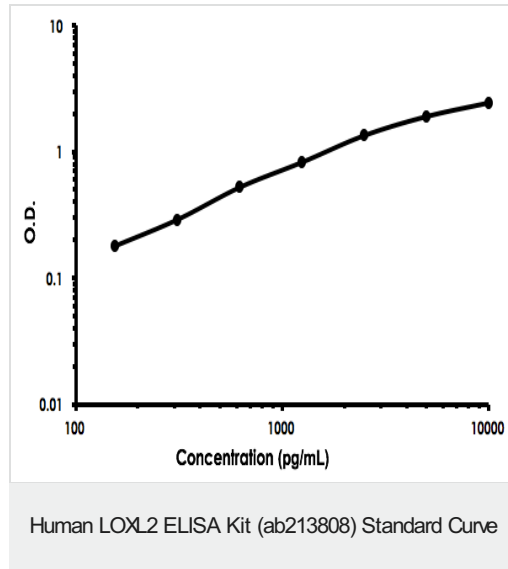
翻訳後修飾

The lysine tyrosylquinone cross-link (LTQ) is generated by condensation of the epsilon-amino group of a lysine with a topaquinone produced by oxidation of tyrosine.

細胞内局在

Secreted > extracellular space.

画像



Human LOXL2 ELISA Kit (ab213808) Standard Curve

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