

DYKDDDDK-Tag Protein ELISA Kit ab285234

2 Images

Overview

Product name	DYKDDDDK-Tag Protein ELISA Kit
Detection method	Colorimetric
Sample type	Cell Lysate
Assay type	Competitive
Sensitivity	>= 5 nM
Range	5 nM - 500 nM
Assay time	1h 30m
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Species independent
Product overview	This DYKDDDDK-Tag Protein ELISA Kit (ab285234, E4700) is based on the competitive ELISA principle and it is an easy, fast and sensitive method to detect the expressed DYKDDDDK-tag proteins. This detection kit offers ready-to-use reagents, and can detect as low as 5 nM of DYKDDDDK-tag protein in approximately 1.5 hrs.
Notes	This product is manufactured by BioVision, an Abcam company and was previously called E4700 DYKDDDDK-Tag Protein ELISA Kit. E4700-100 is the same size as the 96 test size of ab285234.
Platform	Microplate (12 x 8 well strips)

Properties

Storage instructionsStore at -20°C. Please refer to protocols.

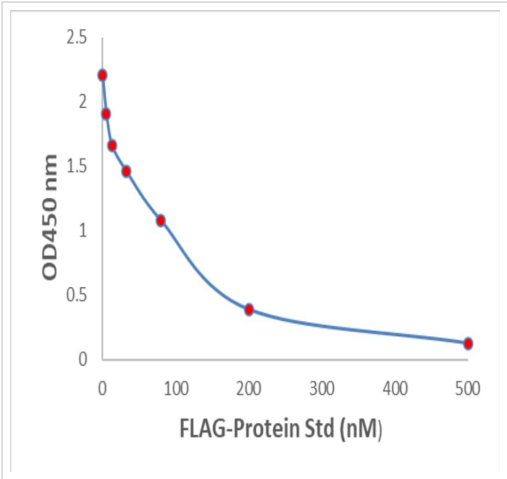
Components	96 tests
10X Wash Buffer II	1 x 50ml
DYKDDDDK Antibody	1 x 7ml
Conjugate Buffer III	1 x 7.5ml
DYKDDDDK Standard	1 vial

Components	96 tests
DYKDDDDK Microplate	1 unit
Goat Anti-Rabbit HRP Conjugate I	1 x 25µl
Microplate Sealing Film	4 units
Stop Solution VIII	1 x 10ml
TMB Substrate I	1 x 10ml

Relevance

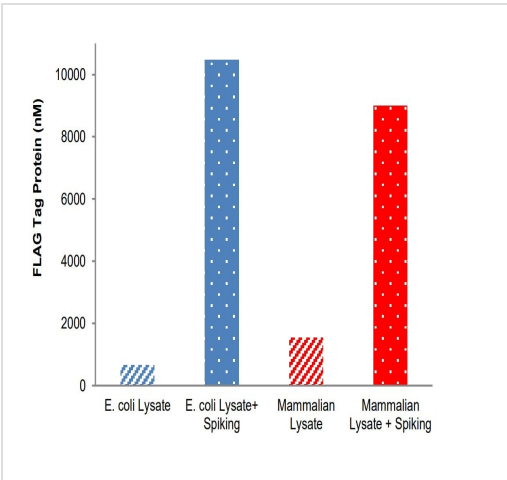
This is a useful tool for the localisation and characterisation of DDDDK tagged proteins (Binds to FLAG® tag sequence).

Images



Standard curve for DYKDDDDK-Tag ELISA Kit (This standard curve is for demonstration only. A standard curve must be run with each assay).

ab285234-DYKDDDDK-Tag-Protein-ELISA-Kit-standard-curve



Spike recovery experiment: E. coli and mammalian cell lysates were spiked with DYKDDDDK-tag protein and showed 88-100% recovery.

ab285234-DYKDDDDK-Tag-Protein-ELISA-Kit-spike-recovery

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