

DNA-Protein Binding Assay Kit (Colorimetric) ab117139

6 References

Overview

Product name DNA-Protein Binding Assay Kit (Colorimetric)

Detection method Colorimetric

Assay time 3h 00m

Species reactivity **Reacts with:** Mouse, Human

Predicted to work with: Mammals 

Product overview

Protein-DNA interactions play a critical role for cellular functions such as signal transduction, gene transcription and epigenetic silencing. Identifying the genetic targets of DNA binding proteins and knowing the mechanism of protein-DNA interaction is important for understanding cellular processes.

DNA-Protein Binding Assay Kit (Colorimetric) (ab117139) allows the user to investigate direct *in vitro* protein-DNA interaction, using dsDNA of choice, within 3 hours. ab117139 is specifically suitable for detection of transcription factor activation in mammalian tissue and cell extracts.

Platform

Microplate reader

Properties

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

Components	96 tests
10X Wash Buffer	1 x 20ml
8 Well Assay Strips (with Frame)	12 units
Antibody Dilution Buffer	1 x 20ml
Assay Binding Buffer	1 x 4ml
Developing Solution	1 x 10ml
Stop Solution	1 x 6ml

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