


DNA-Protein Binding Assay Kit (Fluorometric) ab117140

製品の概要

製品名	DNA-Protein Binding Assay Kit (Fluorometric)
検出方法	Fluorescent
全工程の試験時間	3h 00m
種交差性	交差種: Mouse, Human 交差が予測される動物種: Mammals 

製品の概要

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 (Fluorometric) (ab117140) 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.

試験プラットフォーム

Microplate reader

製品の特長

保存方法

Please refer to protocols.

内容	96 tests
10X Wash Buffer	1 x 20ml
8-Well Assay Strips (with Frame)	12 units
96-Well Assay Plate	1 unit
Antibody Buffer	1 x 6ml
Assay Buffer	1 x 10ml
DTT (500X)	1 x 22µl
Extract Cleaner	1 x 100µl

内容	96 tests
Fluoro Developer	1 x 0.2ml
Oligo Release Solution	1 x 11ml

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