

1 Image

Product name	Mouse DAN ELISA Kit
Detection method	Colorimetric
Precision	Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

**Range** 35 pg/ml - 10000 pg/ml

Recovery	Sample specific recovery
----------	--------------------------

This assay employs an antibody specific for mouse DAN coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures DAN present in the samples. The wells are washed and biotinylated anti-mouse DAN antibody is added. After

washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of DAN bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

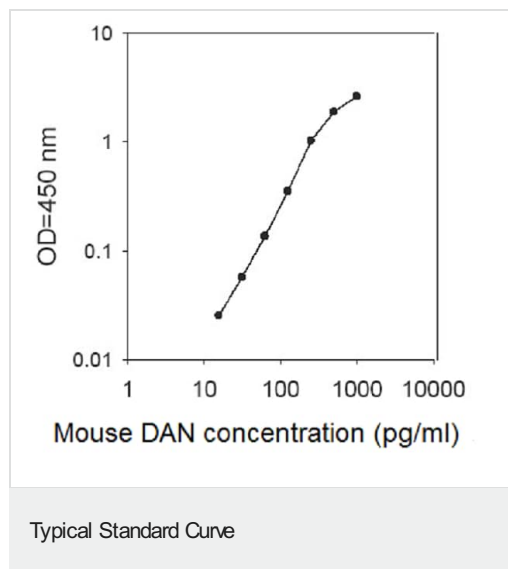
**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
250X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated Mouse DAN detection antibody	2 vials
mouse DAN Standards (lyophilized)	2 vials
Pre-coated mouse DAN Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## Images



Representative Standard Curve using ab193721.

**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