

Standard for TNF alpha Human ELISA Kit ab129028

概述

产品名称	Standard for TNF alpha人ELISA试剂盒
检测方法	Colorimetric
检测类型	Sandwich
实验步骤	Multiple steps standard assay
平台	Microplate

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	1 units
TNF alpha Standard	1 unit

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors