

Product datasheet

Thiol Green Indicator WS ab213601

Overview

Product name Thiol Green Indicator WS

Product overview Ab213601 (Thiol Green Indicator WS) is a component for **ab205811** (GSH/GSSG Ratio Detection Assay Kit II (Fluorometric - Green)). Refer to the kit protocol booklet for instructions of use. Please contact our Scientific Support team for more lot specific information.

Properties

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

Components	1 units
Thiol Green Indicator WS	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.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