

Product datasheet

FirePlex®-HT Imager Setup Kit ab229130

FirePlex®

Overview

Product name	FirePlex®-HT Imager Setup Kit
Detection method	Fluorescent
Assay type	Multiplex
Product overview	<p>FirePlex-HT Imager Setup Kit (ab229130) is designed to identify properly resolving high-content imaging instruments, which can be validated for use with FirePlex-HT® immunoassays:</p> <ul style="list-style-type: none">- Human Cytokines FirePlex®-HT Panel 1 (10-plex) (ab234897)- Human Cytokines FirePlex®-HT Panel 2 (6-plex) (ab234898)- Human Cardiotoxicity FirePlex®-HT Panel 1 (ab252375)- Mouse Cytokines FirePlex®-HT Panel 1 (ab252376)- TGF Beta FirePlex®-HT Panel (ab270739)

The kits contains 384 Well Imaging Plate, plate lid and seal, and HCI Imager Set up Particles

Notes	FirePlex® is a registered trade mark in the United States and is registered as a European Union Trade Mark.
--------------	---

Platform	Multiplex
-----------------	-----------

Properties

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

Components	10 tests
384 Well Imaging Plate	1 unit
384 Well Imaging Plate Lid	1 unit
384 Well Imaging Plate Seal	1 unit
HCI Imager Set up Particles	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