

# HIF1a + PDK1 Hypoxia Response Human Flow Cytometry Kit

## ab126700

5 图像

概述

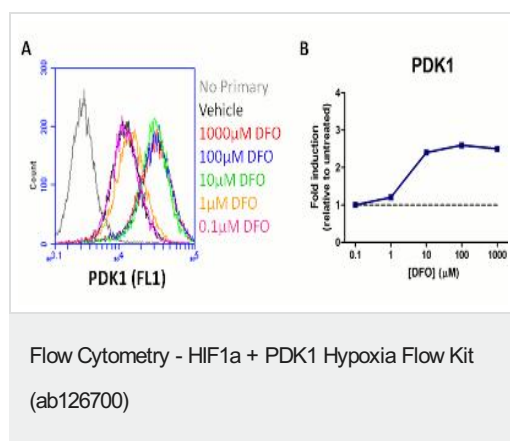
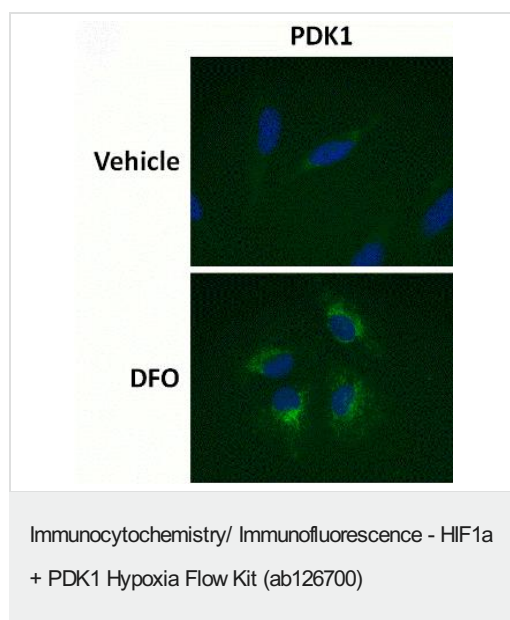
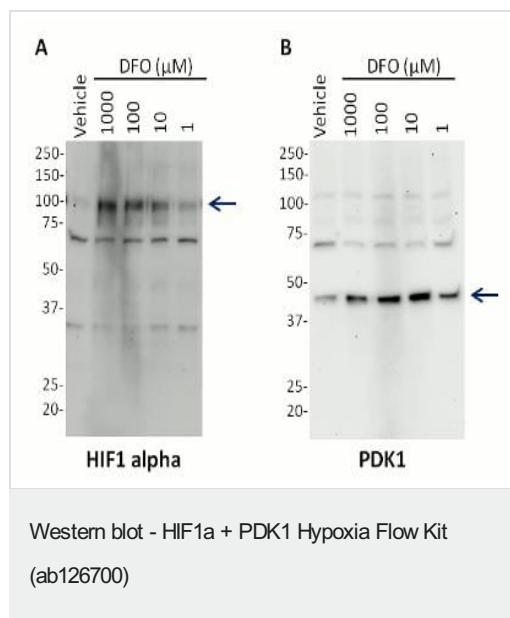
产品名称	HIF1a + PDK1 Hypoxia Response人Flow Cytometry试剂盒
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Sandwich
种属反应性	与反应: Human
产品概述	Hypoxia and the cellular response to hypoxic environment are central topics in studies of metabolism, cancer progression and development and stem cells. A key player is the transcription factor HIF1 alpha (hypoxia inducible factor 1 alpha) which is stabilized at the protein level in response to decreased oxygen tension. HIF1 alpha then promotes transcription of a number of factors that alters cellular physiology. This Hypoxic Response Human Flow Cytometry Kit provides duplexed measurements of the transcription factor HIF1 alpha and the HIF1 alpha responsive protein PDK1.
平台	Flow cytometer

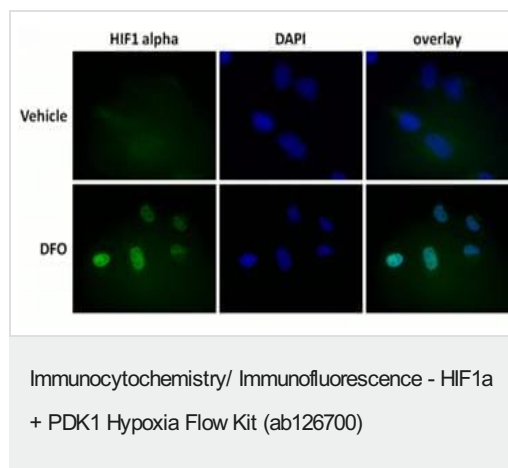
性能

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

组件	1 x 96 tests
100X HIF1A Primary Monoclonal Antibody (Rabbit)	1 x 120µl
100X PDK1 Primary Antibody (Mouse Monoclonal)	1 x 120µl
10X Blocking Buffer	1 x 80ml
10X Phosphate Buffered Saline	1 x 100ml

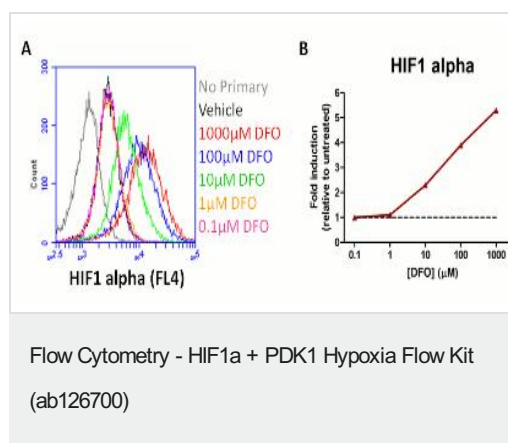
图片





**Figure 3. Antibody specificity demonstrated by**

**immunocytochemistry.** Primary antibodies used in this assay kit were validated by staining HeLa cells +/- treatment with 1mM DFO (24h) and imaged by fluorescent microscopy. HIF1 alpha staining is absent in untreated cells and induced by DFO treatment. HIF1 alpha localizes to the nucleus (as seen by co-localization with the DNA stain DAPI) as expected.



**Figure 1. Sample experiment using ab126700 on HeLa cells treated with a titration of DFO: HIF1 alpha readout.** HeLa cells

were cultured in standard tissue culture plates and treated with a titration of DFO. After 24 hours of DFO exposure, the cells were harvested, fixed and stained as described in the protocol. **(A)** Flow cytometry histogram showing mean fluorescent intensity of HIF1 alpha staining for untreated (Vehicle) and DFO treated samples. In this experiment anti-rabbit-DyLight®650 ([ab96902](#), 1:2000) was used as the secondary antibody and the signal was collected in FL4. **(B)** Plot showing fold induction of HIF1 alpha levels (relative to untreated cells) as a function of DFO concentration (red line). The gray dotted line demarks 1 (the untreated level). DFO concentrations =10µM induce HIF1A protein levels in a dose dependent manner.

**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