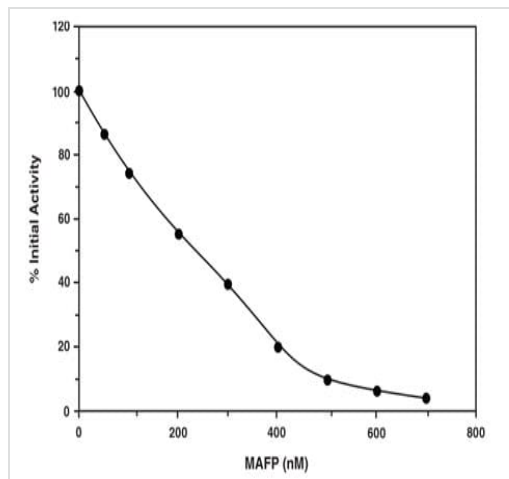


PAF Acetylhydrolase Inhibitor Screening Assay Kit ab133091

1 References 1 图像

概述															
产品名称	PAF Acetylhydrolase抑制剂Screening Assay试剂盒														
检测方法	Colorimetric														
检测类型	Enzyme activity														
产品概述	ab133091 uses 2-thio PAF as a substrate for PAF Acetylhydrolase (PAF-AH). Upon hydrolysis of the acetyl thioester bond at the sn-2 position by PAF Acetylhydrolase, free thiols are detected using 5,5'-dithio-bis-(2-nitrobenzoic acid) (DTNB). The PAF Acetylhydrolase Inhibitor Screening assay includes Human plasma PAF Acetylhydrolase as control and is a time saving tool for screening vast numbers of inhibitors.														
平台	Microplate reader														
性能															
存放说明	Store at -20°C. Please refer to protocols.														
<table><tr><th>组件</th><th>96 tests</th></tr><tr><td>2-thio PAF (substrate)</td><td>2 vials</td></tr><tr><td>96-Well Plate (Colorimetric Assay)</td><td>1 unit</td></tr><tr><td>96-Well Plate Cover</td><td>1 unit</td></tr><tr><td>Human Plasma PAF-AH</td><td>2 vials</td></tr><tr><td>PAF-AH Assay Buffer (10X)</td><td>1 vial</td></tr><tr><td>PAF-AH Assay DTNB</td><td>4 vials</td></tr></table>		组件	96 tests	2-thio PAF (substrate)	2 vials	96-Well Plate (Colorimetric Assay)	1 unit	96-Well Plate Cover	1 unit	Human Plasma PAF-AH	2 vials	PAF-AH Assay Buffer (10X)	1 vial	PAF-AH Assay DTNB	4 vials
组件	96 tests														
2-thio PAF (substrate)	2 vials														
96-Well Plate (Colorimetric Assay)	1 unit														
96-Well Plate Cover	1 unit														
Human Plasma PAF-AH	2 vials														
PAF-AH Assay Buffer (10X)	1 vial														
PAF-AH Assay DTNB	4 vials														
功能	Exhibits lysophospholipase, transacylase, PAF acetylhydrolase and asparaginase activities.														
序列相似性	In the N-terminal section; belongs to the asparaginase 1 family. Contains 5 ANK repeats.														

图片



Inhibition of Human plasma PAF Acetylhydrolase by MAFP (IC_{50} = 250 nM).

Functional Studies - PAF Acetylhydrolase Inhibitor
Screening Assay Kit (ab133091)

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