

Formate Assay Kit ab111748

3 References **1 图像**

概述

产品名称	Formate Assay试剂盒
检测方法	Colorimetric
样品类型	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate
检测类型	Quantitative
检测时间	1h 00m
产品概述	In Abcam's Formate Assay Kit, formate is oxidized to generate a product resulting in color formation ($\lambda = 450 \text{ nm}$) proportional to formate concentration. The kit provides a convenient means for detecting formate in biological samples such as in serum or plasma, cells, culture and fermentation media. There is no need for pretreatment or purification of samples.

Visit our [FAQs page](#) for tips and troubleshooting.

说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K653 Formate Colorimetric Assay Kit. K653-100 is the same size as the 100 test size of ab111748.</p> <p>Formic acid (HCOOH) is the simplest carboxylic acid and occurs naturally, notably in ant and bee venom. It is normally present at low levels in blood and urine (up to $\sim 0.5 \text{ mM}$) but can be present in concentrations as high as 5 mM in acute methanol poisoning. Environmental exposure to formaldehyde can also elevate the blood and urine levels of formate.</p>
----	---

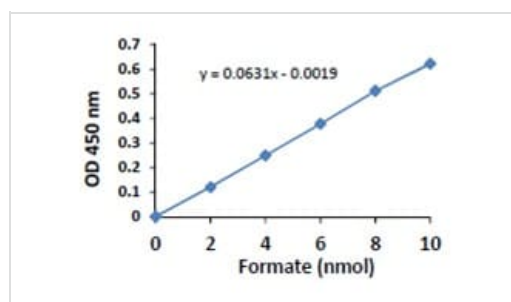
平台	Microplate reader
----	-------------------

性能

存放说明 Store at -20°C . Please refer to protocols.

组件	100 tests
Assay Buffer II	1 x 25ml
Developer Solution III	1 vial
Enzyme Mix IX	1 vial
Formate Standard	1 x 100 μl

图片



Example of standard curve obtained using ab111748

Functional Studies - Formate Assay Kit (ab111748)

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