

Dihydrofolate Reductase Assay Kit (Colorimetric)

ab239705

[1 References](#) [4 Images](#)

Overview

Product name	Dihydrofolate Reductase Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Tissue, Adherent cells, Suspension cells
Product overview	Dihydrofolate Reductase Assay Kit (Colorimetric) (ab239705) is based on the ability of DHFR to catalyze the oxidation of NADPH. The reaction progress is followed by monitoring the decrease in absorbance at 340 nm. Our assay has been optimized to be carried out in a 96-well plate. It is simple, sensitive and can detect as low as 4 mU/ml in a variety of samples.

Notes	This product is manufactured by BioVision, an Abcam company and was previously called K246 Dihydrofolate Reductase Activity Kit (Colorimetric). K246-100 is the same size as the 100 test size of ab239705.
Platform	Microplate reader

Properties

Storage instructions	Please refer to protocols.
----------------------	----------------------------

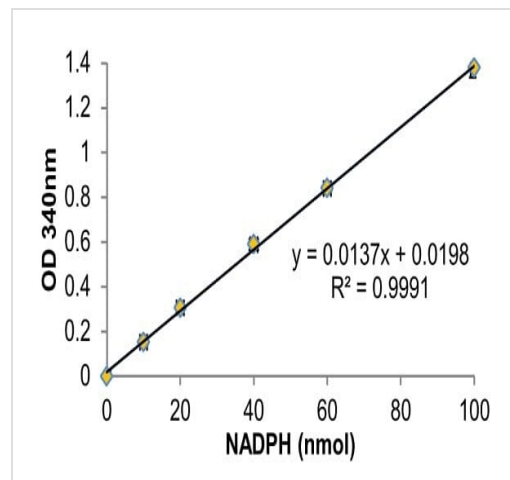
Components	100 tests
DHFR Assay Buffer	1 x 35ml
DHFR Substrate	1 x 450µl
Dihydrofolate Reductase	1 x 10µl
NADPH III	1 vial

Function	Key enzyme in folate metabolism. Catalyzes an essential reaction for de novo glycine and purine synthesis, and for DNA precursor synthesis.
Pathway	Cofactor biosynthesis; tetrahydrofolate biosynthesis; 5,6,7,8-tetrahydrofolate from 7,8-dihydrofolate: step 1/1.

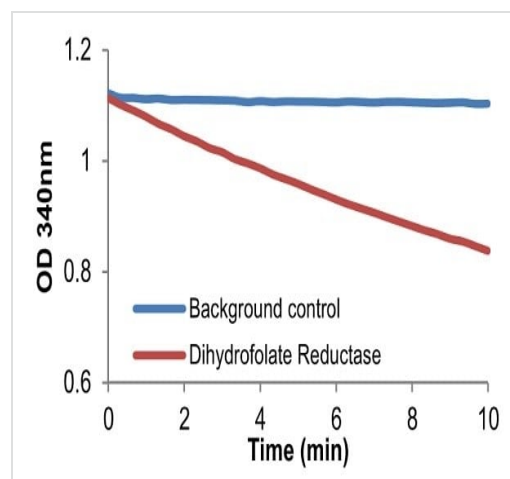
Sequence similarities

Belongs to the dihydrofolate reductase family.
Contains 1 DHFR (dihydrofolate reductase) domain.

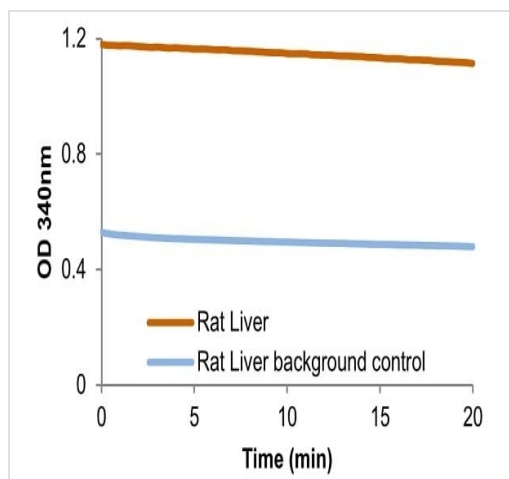
Images



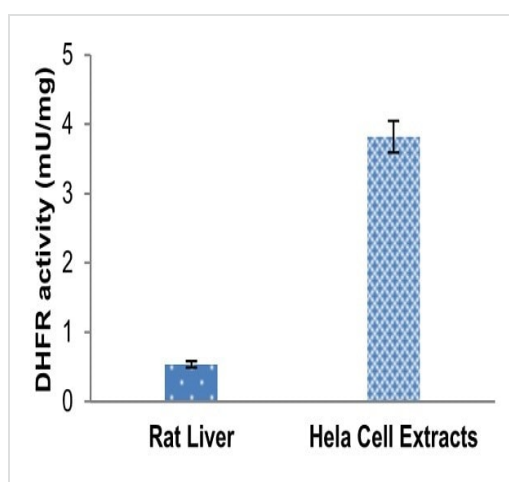
NADPH standard curve



Purified Dihydrofolate Reductase activity



Dihydrofolate Reductase activity in Rat Liver (100 µg)



Dihydrofolate Reductase specific activity was calculated from Rat Liver (100 µg) or HeLa cells (80 µg)

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