

Product datasheet

Sulfatase Activity Assay Kit (Colorimetric) ab204731

[3 Images](#)

Overview

Product name	Sulfatase Activity Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Tissue, Adherent cells, Suspension cells
Assay type	Enzyme activity
Product overview	Sulfatase Activity Assay Kit (ab204731) provides a quick and easy way to determine sulfatase activity. The kit measures the hydrolysis of a sulfate ester to 4-nitrocatechol, which can be detected at OD=515 nm. The kit is suitable for measuring activity of purified enzyme as well as sulfatase from biological samples. The limit of detection is below 1 mU.

Notes

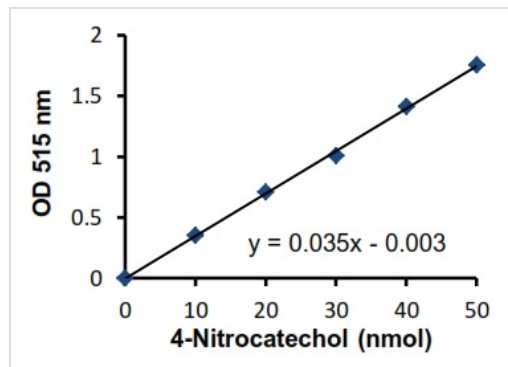
Sulfatases (EC 3.1.6) are enzymes in the esterase class that catalyze the hydrolysis of sulfate esters from a wide range of biological molecules, including steroids, carbohydrates, and proteins. They can be found in intracellular and extracellular spaces and are distributed in a wide range of cells and tissues. Intracellular sulfatases are commonly found localized within the lysosome. Genetic defects in sulfatase can result in certain lysosomal storage disorders and abnormal expression can contribute to certain hormone-dependent cancers, such as breast and prostate cancer.

Platform Microplate reader

Properties

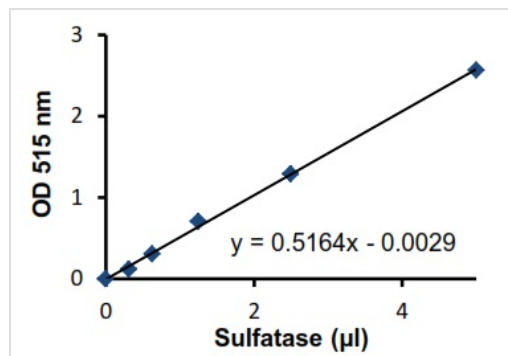
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	100 tests
4-Nitrocatechol Standard (0.5 mM)		1 x 1.5ml
Stop Solution	Red	1 x 10ml
Sulfatase		1 vial
Sulfatase Assay Buffer		1 x 5ml
Sulfatase Substrate		1 x 4ml



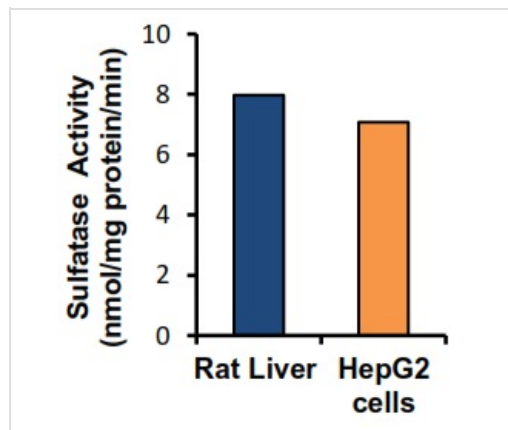
Typical 4-Nitrocatechol Standard calibration curve.

Sulfatase Activity Assay Kit (ab204731)



Enzyme activity measuring different amounts of the provided Sulfatase.

Sulfatase Activity Assay Kit (ab204731)



Measurement of Sulfatase Activity in HepG2 cell lysate (30 µg) and rat liver homogenate (100 µg).

Sulfatase Activity Assay Kit (ab204731)

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