

Arylamine N-acetyltransferase Activity Assay Kit (Fluorometric) ab284533

画像数 3

製品の概要

製品名	Arylamine N-acetyltransferase Activity Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Purified protein
アッセイタイプ	Enzyme activity (quantitative)
ステップ	Multiple steps standard assay
種交差性	交差種: Rat, Human
製品の概要	<p>Arylamine N-acetyltransferase Activity Assay Kit (Fluorometric) (ab284533) (K2031) is a simple, one step plate based assay that measures both NAT1 and NAT2 activity in tissue samples, S9 fractions, recombinant enzyme or purified protein samples.</p> <p>This kit can detect as low as 1 μU in samples.</p> <p>General range: 0-500 pmol/well of acetylated standard.</p>
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K2031 Arylamine N-acetyltransferase Activity Assay Kit (Fluorometric). K2031-100 is the same size as the 100 test size of ab284533.</p>
アプリケーション	適用あり: Enzyme activity assay
試験プラットフォーム	Microplate (12 x 8 well strips)

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Acetylated Standard	1 x 50 μ l
DTT	1 x 100 μ l
NAT Assay Buffer	1 x 25ml

内容	100 tests
NAT Positive Control	1 vial
NAT Substrate I	1 x 200µl
NAT Substrate II	2 vials

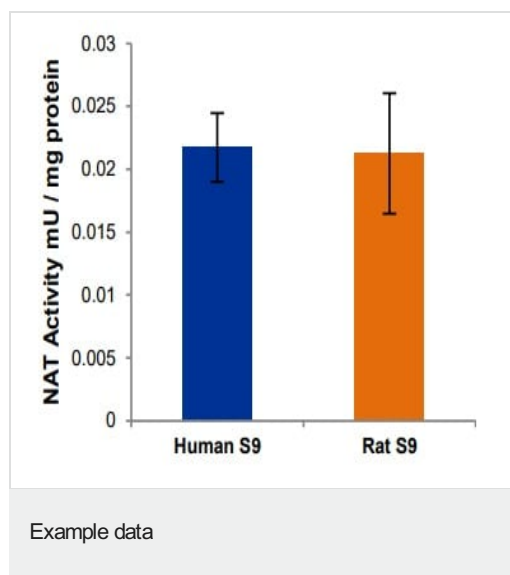
機能	Participates in the detoxification of a plethora of hydrazine and arylamine drugs. Catalyzes the N- or O-acetylation of various arylamine and heterocyclic amine substrates and is able to bioactivate several known carcinogens.
配列類似性	Belongs to the arylamine N-acetyltransferase family.
細胞内局在	Cytoplasm.

アプリケーション

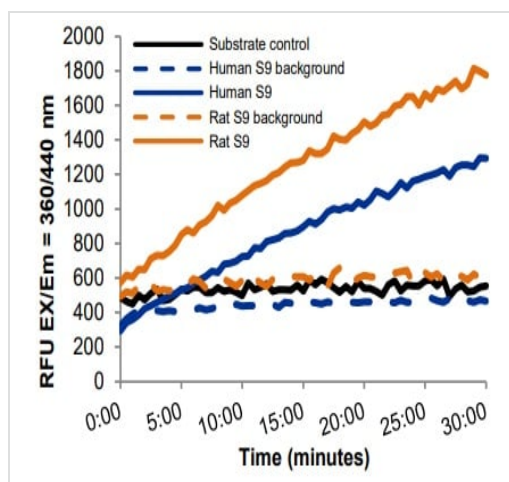
The Abpromise guarantee **Abpromise保証は、**次のテスト済みアプリケーションにおけるab284533の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Enzyme activity assay		Use at an assay dependent concentration.

画像

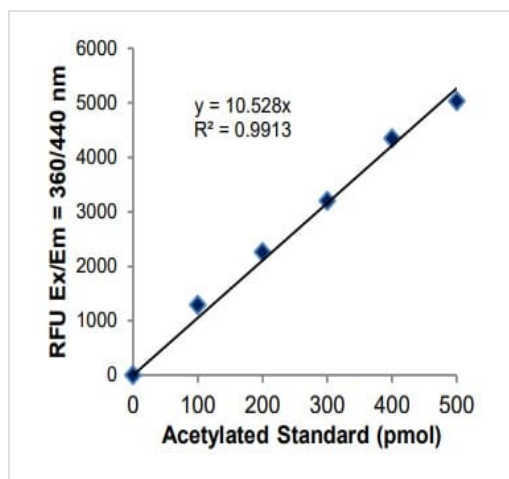


Arylamine N-acetyltransferase specific activity in human and rat liver S9 fractions.



Example data

Enzyme kinetics using Human Liver S9 fraction (160 µg protein/well) and Rat Liver S9 fraction (120 µg protein/well).



Example data

Acetylated Standard Curve. For demonstrative purposes only.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors