

Caspase Family Colorimetric Substrate Kit II Plus ab102488

製品の概要

製品名	Caspase Family Colorimetric Substrate Kit II Plus
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Enzyme activity
製品の概要	<p>Abcam's Caspase Family Colorimetric Substrate Kit II Plus provides ready-to-use colorimetric substrates for members of caspase family proteases. All substrates are provided in liquid ready-to-use form.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K138 Caspase Colorimetric Substrate Set II Plus. K138-9-25 is the same size as the 25 test size of ab102488.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

製品の特性

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

内容	25 tests
2X Reaction Buffer	1 x 20ml
Caspase 6 Substrate, Ac-VEID-PNA	1 x 125µl
Caspase-1 Substrate, Ac-YVAD-pNA	1 x 125µl
Caspase-10 Substrate, Ac-AEVD-PNA	1 x 125µl
Caspase-2 Substrate, Ac-VDVAD-pNA	1 x 125µl
Caspase-3 Substrate, Ac-DEVD-pNA	1 x 125µl
Caspase-4 Substrate, Ac-LEVD-PNA	1 x 125µl

内容	25 tests
Caspase-5 Substrate, Ac-WEHD-pNA	1 x 125µl
Caspase-8 Substrate, Ac-IETD-pNA	1 x 125µl
Caspase-9 Substrate, Ac-LEHD-pNA	1 x 125µl
Cell Lysis Buffer	1 x 100ml
Dilution Buffer	1 x 200ml
DTT	1 x 400µl

関連性 Caspases are members of the cysteine-aspartic acid protease (caspase) family. Sequential activation of caspases plays a central role in the execution-phase of cell apoptosis. Caspases exist as inactive proenzymes which undergo proteolytic processing at conserved aspartic residues to produce 2 subunits, large and small, that dimerize to form the active enzyme.

細胞内局在 Cytoplasmic

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