

Cathepsin G Inhibitor Screening Assay Kit (Colorimetric) ab204693

1 References [画像数 1](#)

製品の概要

製品名	Cathepsin G Inhibitor Screening Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Inhibitor compounds
アッセイタイプ	Enzyme activity
製品の概要	Cathepsin G Inhibitor Screening Assay Kit (Colorimetric) (ab204693) uses the ability of active Cathepsin G to cleave a synthetic pNA (4-Nitroaniline)-based peptide substrate to release pNA, which can be easily quantified using a colorimetric microplate reader. In the presence of a Cathepsin G inhibitor, the cleavage of this substrate is reduced/abolished resulting in decrease or total loss of the pNA absorbance. This simple and high-throughput adaptable assay kit can be used to screen/study/characterize potential inhibitors of Cathepsin G.

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K162 Cathepsin G Inhibitor Screening Kit (Colorimetric). K162-100 is the same size as the 100 test size of ab204693.

Cathepsin G (CTSG, EC 3.4.21.20) is a Serine protease found in azurophilic granules of neutrophilic polymorphonuclear leukocytes. The encoded protease has specificity similar to that of chymotrypsin C, and may participate in the killing and digestion of engulfed pathogens, and in connective tissue remodeling at sites of inflammation.

試験プラットフォーム

Microplate reader

製品の特性

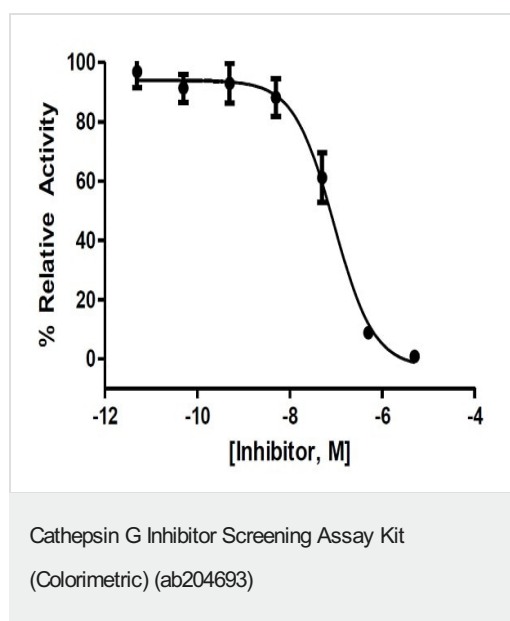
保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Cathepsin G Inhibitor	1 x 20µl
Cathepsin G Substrate	1 x 200µl
CTSG Assay Buffer	1 x 20ml

内容	100 tests
Human Cathepsin G II	1 vial

機能	Serine protease with trypsin- and chymotrypsin-like specificity. Cleaves complement C3. Has antibacterial activity against the Gram-negative bacterium <i>P.aeruginosa</i> , antibacterial activity is inhibited by LPS from <i>P.aeruginosa</i> , Z-Gly-Leu-Phe-CH ₂ Cl and phenylmethylsulfonyl fluoride.
配列類似性	Belongs to the peptidase S1 family. Contains 1 peptidase S1 domain.
細胞内局在	Cell surface.

画像



Inhibition of Cathepsin G activity by CTSG Inhibitor.

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