

Apoptosis Inducer Kit (Ready-to-use) ab102480

5 References

医薬用外劇物

製品の概要

製品名	Apoptosis Inducer Kit (Ready-to-use)
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Direct
製品の概要	Apoptosis Inducer Kit (ab102480) contains 5 ready-to-use chemical reagents that induce apoptosis through different mechanisms. Actinomycin D, an antineoplastic antibiotic, inhibits RNA synthesis. Camptothecin, an inhibitor of nuclear topoisomerase, induces apoptosis in many types of cells. Cycloheximide, an active antibiotic against many yeast and fungi, inhibits protein synthesis. Dexamethasone, an active and highly stable glucocorticoid, probably induces apoptosis by binding and activating the intracellular glucocorticoid receptor. Etoposide, a derivative of podophyllotoxin, inhibits topoisomerase activity. Visit our FAQs page for tips and troubleshooting.
特記事項	This product is manufactured by BioVision, an Abcam company and was previously called K121 Ready-to-use Apoptosis Inducer Set. K121-5 is the same size as the 5 test size of ab102480.

製品の特性

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

内容	5 tests
Actinomycin D 10mM	1 x 50µl
Camptothecin 2mM	1 x 1ml
Cycloheximide 100mM	1 x 1ml
Dexamethasone 10mM	1 x 1ml
Etoposide 100mM	1 x 100µl

関連性 Apoptosis Inducers are chemical reagents that induce apoptosis through various mechanisms.

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