

Apoptosis Inducer Kit (Ready-to-use) ab102480

5 References

概述

产品名称	Apoptosis Inducer试剂盒(Ready-to-use)
样品类型	Adherent cells, Suspension cells
检测类型	Direct
产品概述	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.
说明	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.

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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

相关性 Apoptosis Inducers are chemical reagents that induce apoptosis through various mechanisms.

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.cn/abpromise> or contact our technical team.

Terms and conditions

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