

DNase I Assay Kit (Fluorometric) ab234056

★★★★★ [1 Abreviews](#) [5 References](#) [3 图像](#)

概述

产品名称	DNase I Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Serum, Purified protein
检测类型	Enzyme activity (quantitative)
种属反应性	与反应: Mammals, Other species
产品概述	DNase I Assay Kit (Fluorometric) (ab234056) allows for quantitative evaluation of DNase I activity of purified enzymes and their inhibitors as well as comparative examination of DNase I activity in biological samples. Enzyme activity is detected upon cleavage of a DNA Probe, which yields a fluorescent DNA product measured at Ex/Em = 651/681 nm. The limit of quantification (L.O.Q) is 178 fmoles of DNA probe cleaved per minute per mL.

说明 This product is manufactured by BioVision, an Abcam company and was previously called K429 DNase I Activity Assay Kit (Fluorometric). K429-100 is the same size as the 100 test size of ab234056.

Deoxyribonuclease I (DNase I) is an endonuclease that cleaves DNA phosphodiester bonds yielding 5'-phosphorylated and 3'-hydroxylated oligonucleotides. DNase I targets single-stranded DNA, double-stranded DNA, and chromatin in a non-specific manner. As an important player in cellular waste management, DNase I is normally secreted extracellularly to clear the system from circulating cell-free DNA, foreign DNA from food digestion or potential pathogens, and endogenous chromosomal DNA from apoptotic and necrotic cells. Abnormal DNase I activity occurs in association with a variety of cancers and auto-immune illnesses that exhibit elevated levels of cell-free DNA. Furthermore, DNase I has been therapeutically used in cystic fibrosis patients to degrade DNA and reduce sputum viscosity.

平台 Microplate reader

性能

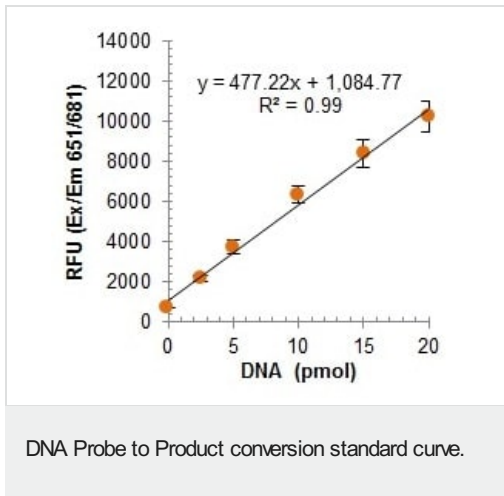
存放说明 Store at -20°C. Please refer to protocols.

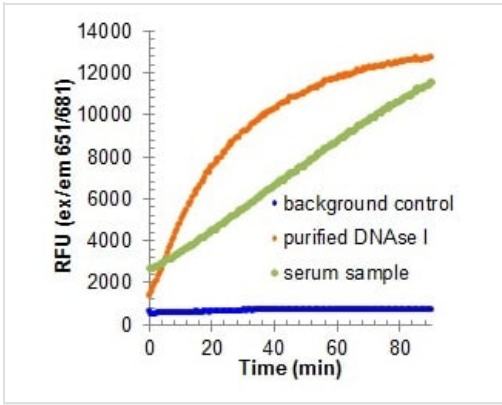
组件	100 tests
10X DNase I Assay Buffer	1 x 1.1ml

组件	100 tests
DNA Probe I	1 vial
DNA Probe Resuspension Buffer	1 x 250µl
DNase I Positive Control	1 vial
Molecular Biology Grade Water	1 x 25ml
Positive Control Resuspension Buffer	1 x 1ml

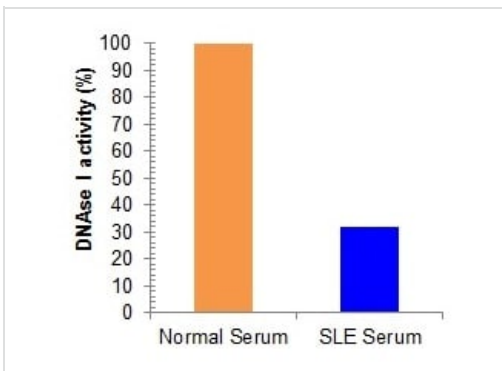
功能	Among other functions, seems to be involved in cell death by apoptosis. Binds specifically to G-actin and blocks actin polymerization.
组织特异性	Principally in tissues of the digestive system. Highest levels found in urine, but also relatively abundant in semen and saliva.
序列相似性	Belongs to the DNase I family.
细胞定位	Secreted. Nucleus envelope. Secretory protein, stored in zymogen granules and found in the nuclear envelope.

图片





Representative activity curve for purified DNase I (orange), serum sample (green), and background control (blue) at 37°C.



Comparative analysis of DNase I activity from 25 µl undiluted single donor normal vs. Systemic Lupus Erythematosus (SLE) patient serum sample.

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