

DNA Isolation Kit - Plasma/Serum ab156893

4 References 1 图像

概述

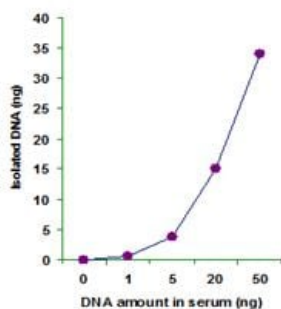
产品名称	DNA分离试剂盒- Plasma/Serum
样品类型	Serum, Plasma, Other biological fluids
范围	0.1 ng - 2000 ng
检测时间	0h 15m
产品概述	<p>DNA Isolation Kit - Plasma/Serum (ab156893) is a complete set of essential components which enables the experimenter to efficiently isolate DNA from plasma, serum, and body fluids. The entire procedure can be completed within 15 minutes and produces far superior results than any competitor kits.</p> <p>ab156893 allows isolation of DNA sizes from 100 bp to 20 kb; DNA quantity from 0.1 ng to 2 µg, optimal at between 1 ng and 1 µg. The isolation uses non-toxic reagents and no phenol chloroform.</p>

性能

存放说明 Please refer to protocols.

组件	100 tests	50 tests
F-Collection Tube	1 x 100 units	1 x 50 units
F-Spin Column	1 x 100 units	1 x 50 units
PS1 (Digestion Solution)	2 x 1.1ml	1 x 1.1ml
PS2 (Digestion Powder)	2 vials	1 vial
PS3 (DNA Isolation Buffer)	2 x 26ml	1 x 26ml
PS4 (DNA Elution Solution)	1 x 2ml	1 x 1ml

图片



Demonstration of DNA Recovery

After treatment with DNA digestion buffer, the DNA is easily recovered in 8-20  $\mu$ l by our specially designed F-Spin Column. DNA is ready for down-stream application.

**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