

Circulating DNA Quantification Kit ab156898

1 References 1 Image

Overview

Product name	Circulating DNA Quantification Kit
Detection method	Fluorescent
Sample type	Serum, Plasma
Assay type	Quantitative
Sensitivity	>= 10 ng/ml
Range	10 ng/ml - 1000 ng/ml
Assay time	0h 30m
Product overview	Circulating DNA Quantification Kit (ab156898) is a complete set of essential components which enables the experimenter to measure and quantify circulating DNA. The entire procedure can be completed within 30 minutes and produces far superior results than any competitor kits. The kit is suitable for quantifying circulating DNA from plasma and serum.
Platform	Microplate reader

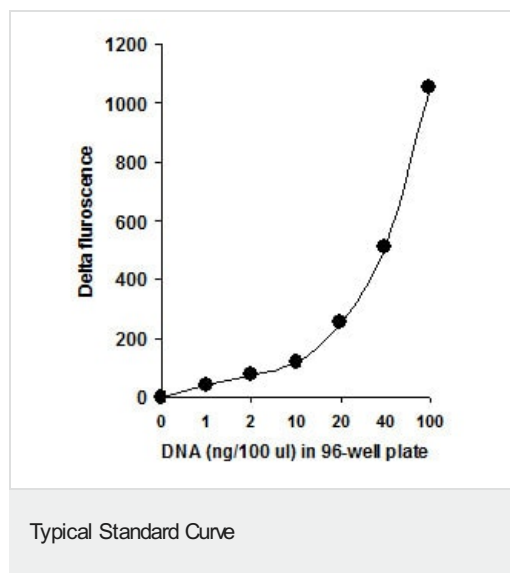
Properties

Storage instructions Please refer to protocols.

Components	48 tests	96 tests
DNQ1 (DNA Digestion Solution)	1 x 1ml	1 x 2ml
DNQ2 (DNA Digestion Powder)	1 vial	2 vials
DNQ3 (DNA Isolation Buffer)	1 x 15ml	1 x 30ml
DNQ4 (DNA Elution Solution)	1 x 1ml	1 x 2ml
DNQ5 (50X DNA Assay Solution)	1 x 0.25ml	1 x 0.5ml
DNQ6 (Assay Dilution Buffer)	1 x 12ml	1 x 24ml
DNQ7 (DNA Standard (10 µg/ml))	1 x 0.1ml	1 x 0.2ml

Components	48 tests	96 tests
F-Collection Tube	1 x 50 units	1 x 100 units
F-Spin Column	1 x 50 units	1 x 100 units

Images



Typical standard curve generated using the Circulating DNA Quantification Kit (ab156898).

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.com/abpromise> or contact our technical team.

Terms and conditions

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