

Cell Fractionation Kit - High Throughput (HT) ab109718

[6 References](#) [3 Images](#)

Overview

Product name	Cell Fractionation Kit - High Throughput (HT)
Assay type	Direct
Assay time	1h 00m
Product overview	Abcam's Cell Fractionation Kit - High Throughput (HT), provides a method and reagents for a simple and rapid preparation of cytosolic, mitochondrial and nuclear fractions from adherent cells.

The Cell Fractionation Kit is designed to produce highly enriched fractions that allow for the monitoring of proteins of interest through various cellular compartments. It is not designed to produce purified cellular fractions.

The kit is especially suitable for, but not limited to, high throughput fractionation of adherent cells in a 96-well plate format. The entire fractionation of 96 samples into the cytosolic, mitochondrial and nuclear fraction is achieved in one hour. The kit is based on sequential detergent extraction of cytosolic, mitochondrial and nuclear proteins without the need for mechanical disruption of cells, and thus fractionates adherent cells into cytosol-containing, mitochondria-containing and nuclei-containing fractions. These fractions are referred throughout the protocol as cytosolic, mitochondrial and nuclear fractions. The kit prepares sufficient sample material for subsequent Western blot analyses, dot blot analyses or for analyses by microplate ELISA assays.

ab109718 is designed to allow the measurement of any proteins which are differentially represented in the cytosol, mitochondria and nuclei, and is particularly applicable to studies of proteins that translocate between these three cellular compartments.

A second version of the kit is also available, [ab109719](#) Cell Fractionation Kit - Standard. This version is designed for fractionating cells grown in batches and is ideal for experiments requiring larger sample amounts.

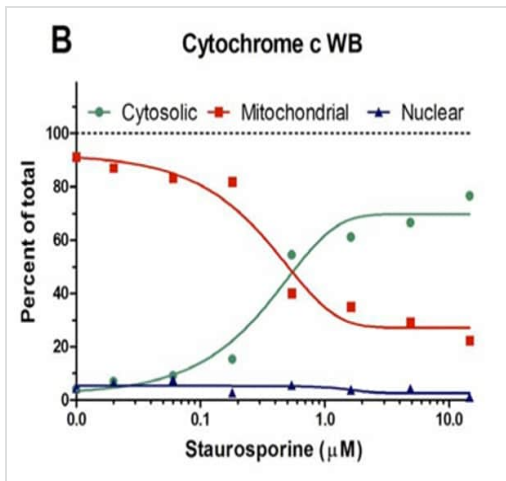
Notes **Storage:** Store Buffer A, Detergent HT II, Detergent HT III and 5X SDS Sample Buffer at -20°C. Store Detergent HT I at -80°C. Store the collection plates and seal films at room temperature.

Properties

Storage instructions Please refer to protocols.

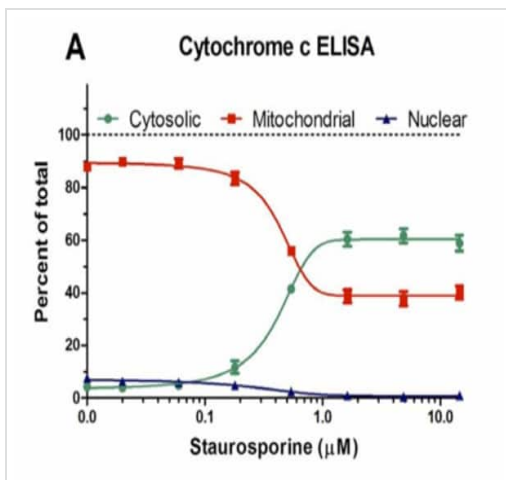
Components	1 x 96 tests
5X SDS Sample Buffer	1 x 1.5ml
96-well collection Plate	3 units
Buffer A	1 x 34ml
Detergent HT I	1 x 7.5µl
Detergent HT II	1 x 35µl
Detergent HT III	1 x 700µl
Plate Seals	3 units

Images



Cell Fractionation Kit - HT (ab109718)

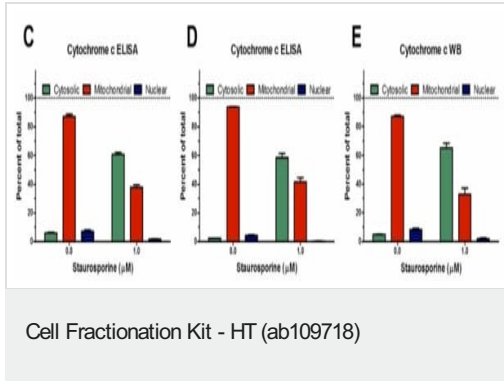
Measurement and distribution of proteins using Western blot, which are differentially represented in the cytosol, mitochondria and nuclei after treatment with varying concentrations of staurosporine.



Cell Fractionation Kit - HT (ab109718)

Measurement and distribution of proteins using an ELISA, which are differentially represented in the cytosol, mitochondria and nuclei after treatment with varying concentrations of staurosporine.

Cell Fractionation Kit - HT (ab109718).



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