

Product datasheet

Aminohippuric acid Assay Kit (Colorimetric) ab204725

1 Image

Overview

Product name	Aminohippuric acid Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Urine, Serum, Plasma
Assay type	Quantitative
Product overview	Aminohippuric acid Assay Kit (ab204725) utilizes a 96-well format in which p-Aminohippuric acid (PAH) reacts with Dimethylaminocinnamaldehyde (DACA) which gives a strongly colored derivative at OD = 550 nm. The amounts of PAH in urine and plasma can easily be quantified with this product.

Notes

p-Aminohippuric acid (PAH) is a derivative of hippuric acid and useful as a diagnostic agent for the measurement of renal plasma flow. About 20-30% is completely filtered by the glomerulus and not reabsorbed by the tubules. The remainder which bypasses the glomerulus and enters the tubules is completely secreted. At low doses, PAH is almost completely removed with one-pass through the kidneys. Hence, the venous concentration of PAH is close to zero and PAH has been used to determine the effective renal plasma flow (eRPF) from the plasma. The venous concentration is usually <10% that of the plasma concentration so eRPF slightly underestimates the actual RPF. This error is generally accepted because of the ease with which PAH infusion allows calculation of eRPF:

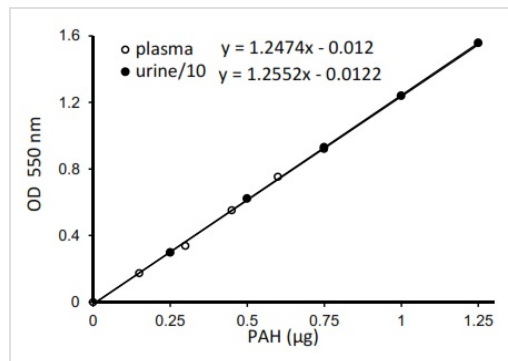
$$eRPF = ([PAH]_U/[PAH]_P) \times \text{Urine Volume}$$

Platform Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
DACA solution	1 x 15ml
PAH Standard (10 mg/ml)	1 x 100µl
TCA (15 %)	1 x 15ml



Typical PAH Standard calibration curve

Aminohippuric acid Assay Kit (Colorimetric)
(ab204725)

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