

## Product datasheet

# Human Follicle Stimulating Hormone ELISA Kit ab108678

### Overview

<b>Product name</b>	Human Follicle Stimulating Hormone ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Serum
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	600 µU/ml
<b>Range</b>	0.001 U/ml - 0.1 U/ml
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	ab108678 Follicle Stimulating Hormone Human ELISA Kit is a immunoenzymatic colorimetric method for quantitative determination of Follicle Stimulating Hormone in serum.

Follicle Stimulating hormone (FSH) is a glycoprotein consisting of two subunits with an approximate molecular mass of 35,500 daltons. The alpha-subunit is similar to other pituitary hormones [luteinizing stimulating hormone (LH), thyroid stimulating hormone (TSH) and chorionic gonadotropin (hCG)] while the beta-subunit is unique. The beta-subunit confers the biological activity to the molecule. Stimulation by gonadotropin-releasing hormone (GnRH) causes release of FSH, as well as LH, from the pituitary and is transported by the blood to their sites of action, the testes or ovary. In men, FSH acts on the Sertoli cells of the testis, stimulating the synthesis of inhibin, which appears to specifically inhibit further FSH secretion, and androgen-binding protein. Thus, it indirectly supports spermatogenesis. In women, FSH acts on the granulosa cells of the ovary, stimulating steroidogenesis. All ovulatory menstrual cycles have a characteristic pattern of FSH, as well as LH, secretion.

<b>Platform</b>	Microplate
-----------------	------------

### Properties

<b>Storage instructions</b>	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	Identifier	1 x 96 tests
10X Wash Solution		1 x 50ml

Components	Identifier	1 x 96 tests
Anti-FSH conjugate		1 x 12ml
Coated Wells	12 strips of 8 wells	1 unit
Foil Cover		1 unit
FSH Control		1 x 1ml
FSH Standard 0 (0 mIU/ml)		1 x 1ml
FSH standard 1 (5 mIU/ml)		1 x 1ml
FSH Standard 2 (10 mIU/ml)		1 x 1ml
FSH Standard 3 (25 mIU/ml)		1 x 1ml
FSH Standard 4 (50 mIU/ml)		1 x 1ml
FSH Standard 5 (100 mIU/ml)		1 x 1ml
Stop Solution		1 x 15ml
Strip holder		1 unit
TMB Substrate Solution		1 x 15ml

**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