

1 References 1 图像

批次内

批次间

特定样本回收率

This product is manufactured by BioVision, an Abcam company and was previously called E4383 BioSim™ Pembrolizumab (Human) ELISA Kit. E4383-100 is the same size as the 96 test size of

ab237652.

平台

Microplate

性能**存放说明**

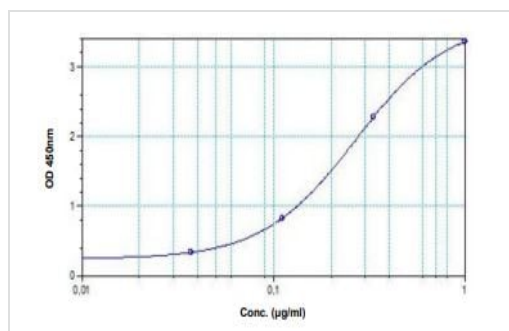
Store at +4°C. Please refer to protocols.

组件	96 tests	96 tests
Assay Buffer	2 x 50ml	2 x 50ml
HRP-conjugate Probe	1 x 12ml	1 x 12ml
Micro ELISA Plate	1 unit	1 unit
Pembrolizumab Standard S1	1 x 0.3ml	1 x 0.3ml
Pembrolizumab Standard S2	1 x 0.3ml	1 x 0.3ml
Pembrolizumab Standard S3	1 x 0.3ml	1 x 0.3ml
Pembrolizumab Standard S4	1 x 0.3ml	1 x 0.3ml
Pembrolizumab Standard S5	1 x 0.3ml	1 x 0.3ml
Pembrolizumab Standard S6	1 x 0.3ml	1 x 0.3ml
Pembrolizumab Standard S7	1 x 0.3ml	1 x 0.3ml
Plate sealers	2 units	2 units
Stop Solution	1 x 12ml	1 x 12ml
TMB substrate	1 x 12ml	1 x 12ml
Wash Buffer (10X)	1 x 50ml	1 x 50ml

相关性

Pembrolizumab is an antibody drug that targets the cell surface receptor programmed cell death protein 1 (PD-1) found on T cells. By preventing the binding of its ligands (PD-L1 and PD-L2), pembrolizumab induces an antitumor immune response. Upregulation of PD-1 ligands is a mechanism for tumors to evade antitumor immune response; when PD-1 binds its ligand, the T cell receives an inhibitory signal leading to T cell anergy and blockade of anti-tumour immune response. Instead of directly targeting tumor tissue to induce tumor cell death, pembrolizumab acts as a checkpoint inhibitor to stimulate immune responses to eliminate cancer cells.

图片



Typical Standard Curve: This standard curve is for demonstration only. A standard curve must be run with each assay.

Example of Pembrolizumab standard curve

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