

OncoSpot™ EGFR T790M FFPE Reference Standard

Catalog number: RM204

Product: EGFR T790M FFPE reference standard

Description: This product can be used as quality control material for optimizing the protocols and evaluating the sensitivity/specificity of assays such as next-generation sequencing (NGS), qPCR and amplification-refractory mutation system (ARMS) PCR. This standard can also be diluted using WT control to generate lower allelic frequencies.

Variant data

AA mutation: T->M

Nucleotide change: ACG ->ATG

Genotype: EGFR (T790M/T790M)

Cosmic ID: COSM6240

Background cell line: HCT116

Allelic Frequency: 100%

Product information

Format: FFPE

Fixation method: 10% neutral buffered formalin solution

Cell density: > 500,000 cells/scroll

Volume/Dimension: 20 µm section

Quantity: 1 scroll

DNA yield: > 400 ng/scroll

Storage: +4 °C

Shelf life: 4 years from the date of manufacture

Quality control: Each lot is tested by sanger sequencing, and H & E staining of 5 µm slide. The DNA extracted from the curl is checked by agarose gel electrophoresis, and Qubit fluorometer. Please refer to the COA for details.



Citation of product: If use of this item results in a publication, please use this information:

OncoSpot™ EGFR T790M FFPE Reference Standard (RM204, GeneCopoeia, Inc., Rockville, MD).

Limited Use License

A limited use license is granted to the Buyer of the Product. The Product shall be used by the Buyer for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products without prior written consent from GeneCopoeia. This Product should be used in accordance with NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Copyright©2019 GeneCopoeia, Inc.

GeneCopoeia, Inc.

9620 Medical Center Drive, #101

Rockville, MD, 20850, USA

Tel: 301-762-0888; Fax: 301-762-3888

Email: support@genecopoeia.com

Web: www.genecopoeia.com