

OncoSpot™ EGFR C797S Genomic DNA Reference Standard

Catalognumber: RM013

Product: EGFR C797S genomic DNA 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.

Variant data

AA mutation: C->S

Nucleotide change: TGC ->AGC

Genotype: EGFR (C797S/C797S)

Cosmic ID: COSM6493937

Background cell line: HCT116

Allelic Frequency: 100%

Product information

Quantity: 1 µg

Concentration: 50 ng/µL

Storage: -20 °C

Shelf life: 3 years from the date of manufacture

Buffer: Tris-EDTA (10 mM Tris-HCl, 1 mM EDTA), pH 8.1

Qualitycontrol: Each lot is tested by sanger sequencing, 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 C797S Genomic DNA Reference Standard, (RM013,
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©2019GeneCopoeia,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