

## **OncoSpot™ KRAS G13D FFPE Reference Standard**

**Catalog number:** RM215

**Product:** KRAS G13D 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:** G->D

**Nucleotide change:** GGC ->GAC

**Genotype:** KRAS (G13D/G13D)

**Cosmic ID:** COSM532

**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™ KRAS G13D FFPE Reference Standard, (RM215, 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](mailto:support@genecopoeia.com)  
Web: [www.genecopoeia.com](http://www.genecopoeia.com)