

Certificate of Analysis

Product name *OncoSpot™ EGFR C797S Genomic DNA Reference Standard*
Product Description *EGFR C797S Genomic DNA Reference Standard*
Zygoty *Homozygote*
Cat. No. *RM013*
Lot. No. *C24WP0101*
Unit Size *1ug (50ng/ul)*

TEST	TEST METHOD	ACCEPTANCE CRITERIA	PASS/FAIL
Genotype	<i>Sanger sequencing of locus specific PCR</i>	<i>Correct base alteration at chromatograph matching the gDNA zygosity</i>	<i>PASS</i>
Quality	<i>Agarose gel electrophoresis</i>	<i>Bright band of high molecular weight</i>	<i>PASS</i>
Quantification	<i>Qubit 4.0 Fluorometer</i>	<i>50±5 ng/μL</i>	<i>PASS</i>

Conclusion **This lot passes the Quality Control specifications.**

Signed by: 

Jiayi Hu

Rockville, MD 20850

USA

Phone: [301-762-0888](tel:301-762-0888) **Fax:** [301-762-3888](tel:301-762-3888)

Website: www.genecopoeia.com

E-Mail: inquiry@genecopoeia.com; support@genecopoeia.com