


Certificate of Analysis

Product name OncoSpot™ EGFR L858R Genomic DNA Reference Standard
Product Description EGFR L858R Genomic DNA Reference Standard
Zygoty Heterozygous
Cat. No. RM003
Lot. No. A22WP0101
Unit Size 1ug (50ng/ul)

TEST	TEST METHOD	ACCEPTANCE CRITERIA	PASS/FAIL
Genotype	Sanger sequencing of locus specific PCR	Correct base alteration at chromatograph matching the gDNA zygosity	PASS
Quality	Agarose gel electrophoresis	Bright band of high molecular weight	PASS
Quantification	Qubit 3.0 Fluorometer	50±5 ng/μL	PASS

Conclusion This lot passes the Quality Control specifications.

Signed by: 
Jiayi Hu

GeneCopoeia, Inc.

Address: 9620 Medical Center Drive, #101
Rockville, MD 20850
USA

Phone: 301-762-0888 **Fax:** 301-762-3888

Website: www.genecopoeia.com

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