

All-in-One™ miRNA Universal Adaptor PCR Primer

Introduction

The **miRNA Universal Adaptor PCR Primer**(Cat.No: QP029) is universal downstream primer for the detection of mature miRNA by adding polyA tail, which is directed at specific olig-dT adaptor design in All-in-One™ miRNA qRT-PCR Detetion Kit 2.0(Cat.No: QP115, QP116), must be used in combination with GeneCopoeia's All-in-One™ miRNA qRT-PCR Detetion Kit 2.0(Cat.No: QP115, QP116), and cannot be applied to other similar miRNA qRT-PCR Detection Kit.

Information

Cat.No	Description	Quantity	Concentration
QP029	Universal Adaptor PCR Primer Tm=64.5 GC content=50%	20 µl	50 µM
	ddH ₂ O(RNase/DNase free)	1 ml	

Storage condition: stored at -20°C, and repeated freezing-thawing is avoided.

Related Products

GeneCopoeia All-in-One™ miRNA qRT-PCR Detetion Kit 2.0(Cat.No: QP115, QP116)
GeneCopoeia All-in-One™ miRNA qPCR Primer (Cat.No: variable).

Procedure

1. Upon receiving the product, the customer shall centrifuge the product at 12000 rpm for 30 sec, so that the liquid flows to the bottom of the tube;
2. This product is the universal downstream primer of specific mature miRNA in the form of 50 µM, which should be diluted by sterilized ddH₂O refer to the PCR reaction system before use;

3. The product should be used in combination with All-in-One™ miRNA qRT-PCR Detection Kit;
4. This product should be used with qPCR reagent containing SYBR Green, not suitable for probe detection.

Application

All-in-One™ miRNA Universal Adaptor PCR Primer is used in combination with **All-in-One™ miRNA qPCR Primer** and **All-in-One™ miRNA qRT-PCR Detection Kit 2.0** to detect the expression of mature miRNA.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2020 GeneCopoeia, Inc.

GeneCopoeia, Inc.
9620 Medical Center Drive, #101
Rockville, MD 20850
USA
Tel: 301-762-0888 Fax: 301-762-3888
Email: inquiry@genecopoeia.com
Web: www.genecopoeia.com