

PRODUCT INFORMATION

pEntranceposon (Kan^R)

#F-766 20 µg

Lot _ **Expiry Date _**

Concentration: 500 ng/µL

Store at -20°C

Description

pEntranceposon (Kan^R) is a high copy plasmid carrying the artificial Mu transposon. The transposon, designated as Entranceposon (Kan^R), has been cloned into the multiple cloning site of the pUC19 vector.

Unit definition

1.0 A₂₆₀ unit dsDNA = 50 µg/mL.

Storage Buffer

10 mM Tris-HCl (pH 8.0), 1 mM EDTA.

Technical support

EMEA: ts.molbio.eu@thermofisher.com

Americas & APAC: ts.molbio@thermofisher.com

Web: www.thermoscientific.com/onebio

PRODUCT USE LIMITATION

This product has been developed and is sold exclusively for research purposes and in vitro use only. This product has not been tested for use in diagnostics or drug development, nor are they suitable for administration to humans or animals. Please refer to www.thermoscientific.com/onebio for Material Safety Data Sheet of the product.

© 2013 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.