



## **REACT<sup>®</sup> 9 Buffer**

**Cat. No. 16316-010**

**Size: 2 × 1 ml**

**Conc.: 10X**

**Store at -20°C or at 4°C.**

For use with Invitrogen restriction endonucleases; detailed instructions are provided on each restriction endonuclease product profile.

Components:

Y90016 REACT 9 Buffer

Buffer Composition (1X concentration):

20 mM Tris acetate (pH 7.9)

10 mM Magnesium acetate

50 mM Potassium acetate

REACT buffers do NOT contain the labile component dithiothreitol (DTT).

REACT buffers do NOT contain bovine serum albumin (BSA). BSA is not necessary in the reaction buffers for Invitrogen restriction endonucleases.

REACT buffer is supplied as the 10X concentrate and should be diluted, 1:10 (1 part REACT buffer + 9 parts other components = 10 parts final reaction volume.)

Doc. Rev.: 052201

This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen Tech-LineSM U.S.A. 800 955 6288

Quality Control:

Endonuclease Assay:

Cleavage by an appropriate restriction endonuclease only at the recognition site at 100-fold excess digest of substrate DNA.

Exonuclease Assay:

Using 1 pmol radiolabeled termini in a 50- $\mu$ l reaction and up to 20 units of an appropriate restriction endonuclease for one hour, the following results were observed:

1 unit of enzyme removed  $\leq$  0.3% label from 5' ends.

1 unit of enzyme removed  $\leq$  0.3% label from 3' ends.

REFER TO THE GIBCO BRL CATALOGUE AND REFERENCE GUIDE  
FOR NOTES ON CONDITIONS WHICH AFFECT ENZYME ACTIVITY.