

Related products

Product	Amount	Cat. no.
Anza™ T4 DNA Ligase Master Mix	50 reactions	IVGN210-4
Anza™ Alkaline Phosphatase Kit	500 reactions	IVGN220-4
Anza™ T4 PNK Kit	500 reactions	IVGN230-4
Anza™ DNA Blunting Kit	100 reactions	IVGN240-4
Anza™ DNA End Repair Kit	20 reactions	IVGN250-4
ReadyPouch™ 1% Agarose Gels	10 pouches	A24574
BlueJuice™ Gel Loading Buffer (10X)	3 x 1 mL	10816-015
One Shot™ TOP10 Chemically Competent <i>E. Coli</i>	20 reactions	C4040-03

To order additional Anza™ Restriction Enzymes, go to
thermofisher.com/Anza

Important licensing information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

©2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Anza™ 73 BclI (10 U/μL)

Cat. No.	Size	Lot no.	Exp. Date
IVGN073-6	2500 units		

Publication No. MAN0014200 Rev. A.0

BclI restriction
site



Components	IVGN073-6
Anza™ 73 BclI enzyme	250 μL
Anza™ 10X Buffer	1000 μL
Anza™ 10X Red Buffer	1000 μL

Storage

Store at -20°C.



For support email thermofisher.com/techresources

thermofisher.com/lifescience

21 June 2015

ThermoFisher
SCIENTIFIC

For research use only. Not for use in diagnostic procedures.

Protocol

1. Prepare a reaction mix by adding reagents in the order indicated in Table 1.
2. Incubate at 37°C for 15 minutes.

Table 1

Reagent	1 enzyme reaction	2 enzyme reaction	3 enzyme reaction
Nuclease-Free Water	As required to make up final reaction volume		
Anza™ 10X Buffer or Anza 10X Red Buffer	2 µL	2 µL	3 µL
DNA	0.2–1 µg	0.2–1 µg	0.2–1 µg
Anza™ restriction enzyme 1	1 µL	1 µL	1 µL
Anza™ restriction enzyme 2	—	1 µL	1 µL
Anza™ restriction enzyme 3	—	—	1 µL
Final reaction volume	20 µL	20 µL	30 µL

Volumes can be scaled up linearly to 5X.

Product description

Invitrogen™ Anza™ Restriction Enzymes are part of a complete cloning system centered around the proprietary Anza™ Buffer.

Anza™ restriction enzymes require no more than 15 minutes for complete digestion of any DNA substrate, but also permit digestion up to 16 hours without deleterious effects.

All Anza™ restriction enzymes and Anza™ modifying cloning enzymes are fully functional in Anza™ Buffer.

Anza™ 10X Buffer

Use Anza™ 10X Buffer for applications that require analysis by fluorescence excitation. Add $\frac{1}{10}$ volume of 10X Loading Buffer to samples digested in Anza™ Buffer before performing electrophoresis.

Anza™ 10X Red Buffer

Anza™ 10X Red Buffer includes a density reagent containing red and yellow tracking dyes.

Load samples digested in Anza™ Red Buffer directly onto a gel for electrophoresis.