

# *C. elegans* transcription factors

Cat. #OCE4819, OCE4821, OCE4818, OCE4820

## Product description

The *C. elegans* Transcription Factor Collection is a genomic resource for the discovery of transcription factor–promoter interactions by means of yeast one-hybrid (Y1H) screens (Vermeissen *et al.*, 2007). Each prey construct contains one of 755 plasmid-encoded *C. elegans* transcription factor ORFs fused with the GAL4p activation domain. The expressed fusion-protein product will induce a GAL4-dependent reporter when the transcription factor domain binds to a promoter bait immediately upstream of the reporter (Deplancke *et al.*, 2006). Each construct is available in yeast (Cat #OCE4820 and OCE4821) for mating with bait strains or in an *E. coli* host (Cat #OCE4818 and OCE4819) for transformation into bait strains.

## Plasmid

pDEST<sub>AD</sub>

## Yeast strain

Y1Ha001 (MAT $\alpha$ ) is derived from Y187.

## Genotype

MAT $\alpha$  ura3-52, his3-200, ade2-101, trp1-901, leu2-3, 112, met-, gal4 $\Delta$ , gal80 $\Delta$ , URA3::GAL1<sub>UAS</sub>-GAL1<sub>TATA</sub>-LacZ  
(see Vermeissen *et al.*, 2007 Supplementary Methods for details)

## Product applications

The *C. elegans* Transcription Factor Collection was created to enable the discovery of transcription factor–promoter interactions by means of yeast one-hybrid (Y1H) assays (Vermeissen *et al.*, 2007). Bait reporter strains are not included in this collection, but may be constructed from any of the cloned promoters in the *C. elegans* Promoter Collection (Cat #PCE1181) by recombinational cloning. See Deplancke *et al.*, (2007) for details.

Alternatively, the *C. elegans* Transcription Factor Collection can be used as prey strains in yeast two-hybrid (Y2H) assays for the purpose of discovering homodimer and heterodimer transcription factor complexes. Appropriate bait strains may be constructed from the *C. elegans* ORF Collection (Cat #OCE4518) by recombinational cloning.

## Protocols

1. For information on recombinational cloning using the Gateway™ Cloning System, please visit the Gateway™ Technology webpage on Invitrogen™'s website.
2. For information on constructing promoter bait strains, please see the reference Deplancke *et al.*, 2004.
3. For information on performing library screens, matrix assays or pooled assays please see the reference Vermeissen *et al.*, 2007.

## Protocols

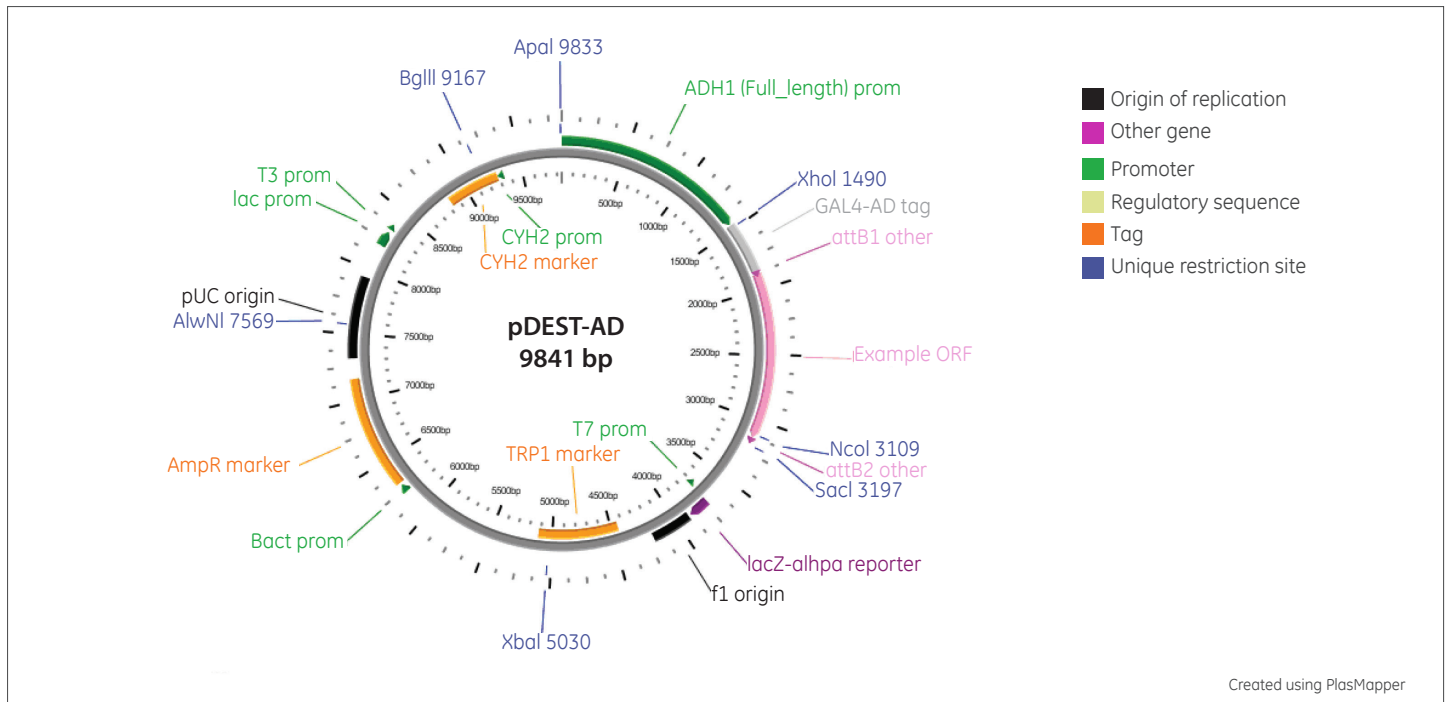


Figure 1. Vector map of pDEST-AD.

## References and suggested reading

1. Deplancke B *et al.*, (2004). A Gateway-compatible yeast one-hybrid system. *Genome Research* **14**, 2093–2101.
2. Deplancke B *et al.*, (2006). A gene-centered *C. elegans* protein-DNA interaction network. *Cell* **125**, 1193–1205.
3. Vermeirssen V *et al.*, (2007). Matrix and Steiner-triple-system smart pooling assays for high-performance transcription regulatory network mapping. *Nature Methods* **4**, 659–664.
4. Walhout *et al.*, (2000). GATEWAY recombinational cloning: application to the cloning of large numbers of open reading frames or ORFeomes. *Methods Enzymol.* **328**, 575–592.

### If you have any questions, contact

**t** +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

**f** + 44 (0)1223 655 581

**w** [horizondiscovery.com/contact-us](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://dharmacon.horizondiscovery.com/service-and-support)

**Horizon Discovery**, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

Gateway and Invitrogen are trademarks of Thermo Fisher Scientific, Inc. ©2018 Horizon Discovery Group Company—All rights reserved. All trademarks are the property of Horizon Discovery Company unless otherwise specified. First published June 2015. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

**horizon**<sup>™</sup>  
INSPIRED CELL SOLUTIONS