

C. elegans transcription factors

Table 1. Applicable catalog numbers

Cat. #	Description
OCE4818	<i>C. elegans</i> Transcription Factor Clone— <i>E. coli</i>
OCE4819	<i>C. elegans</i> Transcription Factor Collection— <i>E. coli</i>
OCE4820	<i>C. elegans</i> Transcription Factor Strain—Yeast
OCE4821	<i>C. elegans</i> Transcription Factor Collection—Yeast

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_{AD}

Yeast strain

Y1Ha001 (MAT α) is derived from Y187.

Genotype

MAT α *ura3-52*, *his3-200*, *ade2-101*, *trp1-901*, *leu2-3*, 112, *met-*, *gal4 Δ* , *gal80 Δ* , URA3::GAL1UAS-GAL1TATA-LacZ)
(see Vermeissen *et al*, 2007 Supplementary Methods for details)

Quality control

All cultures are checked for growth prior to shipment.

Shipping and storage

Individual *E. coli* constructs are shipped as bacterial cultures of *E. coli* in 2x LB broth (low salt)* medium with 8% glycerol, 100 μ g/mL ampicillin. Individual yeast constructs are shipped as yeast cultures of Y1Ha001 in SD(-) tryptophan broth with 15% glycerol. Individual constructs are shipped at room temperature. Collections are shipped in 96-well plate format on dry ice. Individual constructs and collections should be stored at -80°C .

To allow any CO₂ that may have dissolved into the medium from the dry ice in shipping to dissipate, please store constructs at -80°C for at least 48 hours before thawing.

*1x LB medium can be used instead of 2x LB broth medium.

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 **or** dharmacon.horizondiscovery.com/service-and-support

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

©2018 Horizon Discovery Group Company—All rights reserved. All trademarks are the property of Horizon Discovery Company unless otherwise specified. First published June 2014. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.