

# YKO parental strains

Cat. #YSC1048, YSC1049, YSC1050, YSC1061, YSC1062

A nearly complete set of yeast open reading frame (ORF) knock-outs has been produced by the *Saccharomyces* Genome Deletion Project (SGDP) using five Yeast Knockout (YKO) Parental Strains derived from *Saccharomyces cerevisiae* S288C<sup>1</sup>. A PCR-based strategy was used to replace each ORF with a KanMX cassette containing unique tags, “barcodes”, for each deletion. Four different mutant collections have been generated: haploids of mating types MATa and MATalpha, homozygous diploids for non-essential genes, and heterozygous diploids, which contain the essential and non essential ORFs.

## Storage

- 4 °C for up to one week
- –80 °C indefinitely

## 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

©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.

## Product description

Culture of *S. cerevisiae* in YPD broth with an inert growth indicator + 15% glycerol.

BY4730	MATa leu2Δ0 met15Δ0 ura3Δ0
BY4739	MATa leu2Δ0 lys2Δ0 ura3Δ0
BY4741	MATa his3Δ1 leu2Δ0 met15Δ0 ura3Δ0
BY4742	MATa his3Δ1 leu2Δ0 lys2Δ0 ura3Δ0
BY4743	4741/4742

Homozygous diploids are in the BY4743 background unless 4730/4739 is indicated.

## Useful websites and references

*Saccharomyces* Genome Deletion Project, click [here](#).

1. Winzeler, E.A., *et al.*, Functional characterization of the *Saccharomyces cerevisiae* Genome by gene deletion and parallel analysis, *Science*, **285**, 901-906, 1999.