

## Ubiquitin Conjugating Enzymes (E2) sampler pack

**Product Number:** BML-UW8975

The sampler pack contains 10µg quantities of the following enzymes:

Product	Details	Cat. No.
Conjugating enzyme UbcH1	His-tagged (h, rec)	BML-UW9020
Conjugating enzyme UbcH2	His-tagged (h, rec)	BML-UW9025
Conjugating enzyme UbcH3	His-tagged (h, rec)	BML-UW8730
Conjugating enzyme UbcH5a	His-tagged (h, rec)	BML-UW9050
Conjugating enzyme [ <sup>85</sup> C]Ubc H5a	His-tagged [ <sup>85</sup> C mutant] (h, rec)	BML-UW9055
Conjugating enzyme UbcH5b	His-tagged (h, rec)	BML-UW9060
Conjugating enzyme [ <sup>85</sup> C]UbcH5b	His-tagged [ <sup>85</sup> C mutant] (h, rec)	BML-UW9065
Conjugating enzyme UbcH5c	His-tagged (h, rec)	BML-UW9070
Conjugating enzyme [ <sup>85</sup> C]UbcH5c	His-tagged [ <sup>85</sup> C mutant] (h, rec)	BML-UW9075
Conjugating enzyme UbcH6	His-tagged (h, rec)	BML-UW8710
Conjugating enzyme UbcH7	His-tagged (h, rec)	BML-UW9080
Conjugating enzyme UbcH8	His-tagged (h, rec)	BML-UW9135
Conjugating enzyme UbcH9	un-tagged (h, rec)	BML-UW9320
Conjugating enzyme UbcH10	un-tagged (h, rec)	BML-UW0960
Conjugating enzyme UbcH12	His-tagged (h, rec)	BML-UW9145
Conjugating enzyme UbcH13/Mms2	His-tagged (h, rec)	BML-UW9565