

PRODUCT INFORMATION SHEET

Product name: Dynal SSP *AllSet*⁺™ Negative Control
Product number: 990.03
Volume: 24 tests
Kit Batch number: RK0792
Lot number: 343-x
Expiry date: 2007-04
Storage: 2 - 8°C
Manufactured by: Dynal Biotech Ltd., U.K.

The *Dynal AllSet*⁺™ *Negative Control* test is intended to be run in parallel with Dynal SSP typings to check for contamination that may be present in DNA diluent, buffers or Taq Polymerase. The tube contains primer pairs that will detect contamination with PCR amplified fragments derived from Dynal SSP products amplified via the internal control primers. **Contamination with human genomic DNA will also be detected.**

Each *Dynal AllSet*⁺™ *Negative Control* kit contains 3 PCR strips where each well in the PCR strip contains an identical pre-aliquoted dried SSP primer solution. One tube should be cut off and used per test.

Instructions for Use of the *Dynal AllSet*⁺™ *Negative Control* kit:

Mix 2µl of the diluent normally used to dilute your sample DNA (**minus DNA**) with 8µl of the Dynal Mastermix/DynaMix solution and 0,1µl of Taq polymerase and then add to the dried pre-aliquoted negative control. Where there is **no** contamination, **no** detectable PCR fragment will be seen when run under the same PCR conditions as Dynal SSP sets.

Detail of the SSP protocol is available in the IFU supplied with *Dynal AllSet*⁺™ SSP kits.

In the case of contamination, these contamination control primers will give rise to a 280, 429 or a 840 base pair fragment if the contamination originates from a Dynal SSP kit.

Please note that presence of any fragment from the contamination control indicates contamination.

This product may not detect contamination from other PCR based tissue typing kits.

The *Dynal AllSet*⁺™ *Negative Control* primers amplify a segment of the human growth hormone and the HLA DRA gene.

5' primer GCCTTCCCAACCATTCCTTA
GAGGTAAGTGTGCTCACGAACAGC

3' primers TCACGGATTTCTGTTGTGTTTC
GAGAAAGGCCTGGAGGATTC
GGTCCATACCCCAGTGCTTGAGAAG

