

901-0005 Protocol

HRP Conjugate Stabilizer/Diluent Protocol

Introduction:

The performance of HRP conjugates diminishes over time. There are a number of factors that contribute to loss of performance including:

1. Inactivation of HRP
2. Instability of the bond which links HRP to the antibody
3. Microbial attack
4. Denaturation of the antibody

Loss of performance accelerates with increasing temperature and with increasing dilution of conjugate. For these reasons HRP conjugates are often kept in a concentrated form at 4 degrees Celsius. LifeXtend(TM) conjugate stabilizer/diluent is a proprietary multi-component reagent system that protects antibody-HRP conjugates from all of the negative factors noted above, thus ensuring the best possible performance in experiments performed at room temperature. Moreover, the ability to store HRP conjugates at working concentration eliminates waste and improves consistency from experiment to experiment. In addition to its proprietary enzyme stabilizers, LifeXtend(TM) contains antimicrobial agents as well as blockers to enhance signal to background in immunoassay applications.

Storage of LifeXtend(TM):

LifeXtend(TM) HRP conjugate stabilizer/diluent is shipped at ambient temperature and should be stored at 4 degrees Celsius.

Related products:

Lightning-Link HRP conjugation kits - the world's easiest conjugation system, requiring just thirty seconds hands-on time to set up antibody conjugation reactions.

701-0000 LL-HRP conjugation kit, 3 reactions, each up to 400ug of Ab

701-0002 LL-HRP conjugation kit, 1 reaction up to 4mg of Ab Other pack sizes also available.