

Protocol for ChIP-IT® Fixation Buffer

Catalog No.: 53038

Name: ChIP-IT® Fixation Buffer

Format: 2 x 1.5 ml

Store at: 4°C

The following protocol is designed to be used with Active Motif's ChIP-IT® Fixation Buffer for formaldehyde fixation of chromatin for use in chromatin immunoprecipitation (ChIP). This specially formulated buffer can be applied to any chromatin preparation method to provide consistency to the chromatin preparation, resulting in better ChIP results.

Preparation of Complete Fixation Solution

1. Complete Fixation Solution should be prepared fresh before each experiment. Fixation Solution is used at 1:10 growth medium volume. Follow the recommendations in the table below for preparing Complete Fixation Solution based on the growth medium volume to be used. Volumes include excess for pipetting. Adjust volumes as needed for growth conditions.

Reagents	Growth Medium Volume					
	1 ml medium	2 ml medium	3 ml medium	5 ml medium	10 ml medium	20 ml medium
Sterile water	69 µl	138 µl	207 µl	345 µl	785 µl	1.57 ml
ChIP-IT Fixation Buffer	8 µl	16 µl	24 µl	40 µl	90 µl	180 µl
37% Formaldehyde	33 µl	66 µl	99 µl	165 µl	375 µl	750 µl
Total Volume	110 µl	220 µl	330 µl	550 µl	1.25 ml	2.5 ml

Note: Use appropriate precautions when working with formaldehyde (*i.e.* safety glasses, gloves and lab coat).

2. Vortex to mix solution.
3. Use 1/10 growth medium volume per plate (*i.e.* 2 ml Complete Fixation Solution per 20 ml growth medium).
4. Shake gently at room temperature for 15 minutes.
5. Use the conditions of your preferred chromatin preparation protocol to stop the fixation reaction and proceed with chromatin preparation.