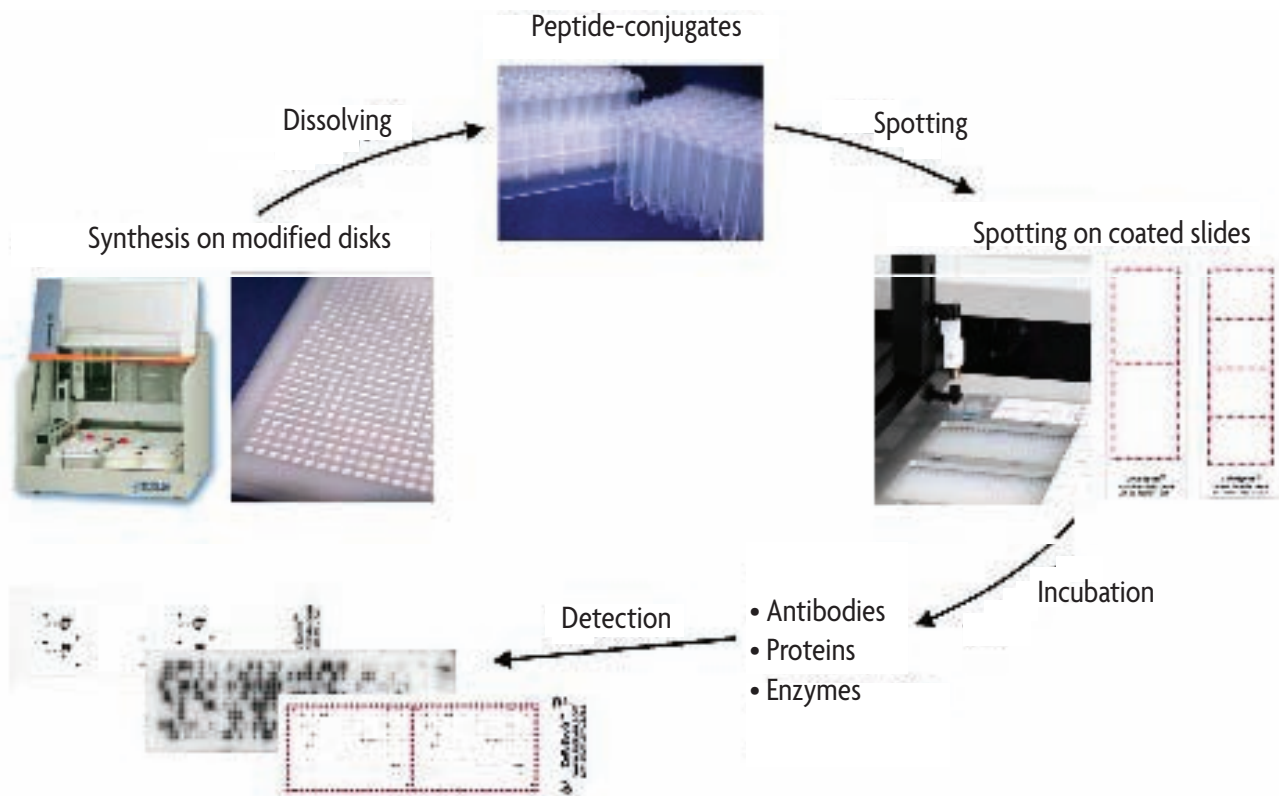


## MODified™ Histone Peptide Arrays

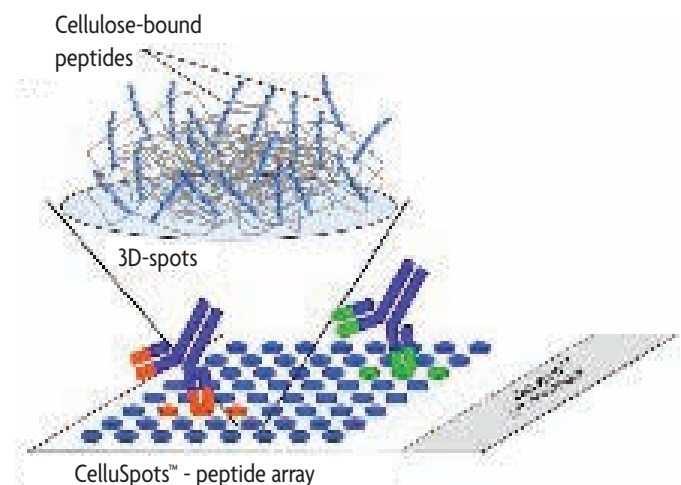
MODified™ Histone Peptide arrays\* are arrays of peptide-cellulose-conjugates spotted on planar surfaces such as coated glass slides. Peptides are synthesized on a modified cellulose support which can be dissolved after synthesis. The solutions of individual peptides covalently linked to macromolecular cellulose are then spotted onto white coated microscope slides. After evaporation of the solvent a three-dimensional layer is formed which is not dissolved in aqueous reagents used for standard assays. The peptide loading in the three dimensional layer of cellulose-peptide-conjugates exceeds that of a monolayer by a factor of 100. This high peptide density in the spots is advantageous to identify even protein-interaction sites with low binding constants

### MODified Histone Peptide Array production:



### Specifications:

- Standard microscope slides (26x76 mm, white coating)
- Up to 768 spots per slide (384 peptide-conjugate spots printed in duplicate)
- Spot-to-spot distance 1.2 mm
- Peptides are covalently bound to cellulose via C-terminus
- Arrays contain control peptides and location marks
- Detection methods include chemiluminescence, chromogenic or autoradiography



\*CelluSpots™ arrays are manufactured under license by INTAVIS Bioanalytical Instruments AG and sold through Active Motif as MODified™ Histone Peptide Arrays.