

abcam

# Gelatin zymography protocol

Detect MMP activity in conditioned media.

Gelatin zymography is a method to detect activity of gelatinase enzymes, such as the matrix metalloproteinases (MMPs) MMP-2 and MMP-9. Active gelatinases digest gelatin embedded in a polyacrylamide gel. After Coomassie staining, areas of degradation are visible as clear bands against a darkly stained background.

This protocol is optimized for detecting secreted MMP-9 and MMP-2 activity in conditioned media.

## Preparation of conditioned media

1. Plate cells cultured in fetal bovine serum (FBS) in a six-well plate (2 mL/well) or in a 75 cm<sup>2</sup> flask (10 mL)
2. At 70–80% confluency, remove FBS media, wash cells twice with RBS-free media and continue to grow cells in FBS-free media.

*The duration of growth in FBS-free media must be optimized for the cell line: for example, 231G and 468 breast cancer cells require a 40–44 h growth period before collection of conditioned media.*

3. Collect conditioned media and centrifuge or filter to eliminate dead cells
4. Concentrate conditioned media 10X

## Gelatin zymography

### Running the gel

1. Adjust conditioned media in all samples to the same protein concentration.

*For each sample, test one aliquot at a low protein concentration (5 µg/mL) and one at a high protein concentration (15 µg/mL).*

2. Add 5x non-reducing sample buffer to your samples
3. Prepare a 7.5% acrylamide gel containing gelatin. Use a 1 mm thickness plate.

Separating gel (7.5 % acrylamide)		Stacking gel	
1.5 M Tris pH 8.8	2 mL	0.5 M Tris pH 6.8	1.25 mL
30% acrylamide	2 mL	30% acrylamide	0.670 mL
H <sub>2</sub> O	2 mL	H <sub>2</sub> O	3.075 mL
Gelatin (4 mg/mL)	2 mL	10 % SDS	50 µL
10% SDS	80 µL	10 % APS	50 µL
10% APS	80 µL	TEMED	10 µL
TEMED	10 µL		

4. Load sample to each well; typically, 10 µL protein per well is suitable. Include a protein molecular weight marker in one well. Run the gel at 150 V until good band separation is achieved.

### Gel washing and staining

1. Wash the gel 2 x 30 min with washing buffer.

*Soaking the gel in washing buffer removes SDS from the gel.*

2. Rinse for 5–10 min in incubation buffer at 37°C with agitation.
3. Replace with fresh incubation buffer and incubate for 24 h at 37°C.

*The incubation buffer contains cofactors necessary for the gelatinase reaction to occur.*

4. Stain the gel with staining solution for 30 min to 1 h. Rinse with H<sub>2</sub>O.
5. Incubate with destaining solution until bands can clearly be seen.

*Areas of enzyme activity appear as white bands against a dark blue background.*

## Buffers

### 5X non-reducing sample buffer

Final concentration	For 250 mL
4% SDS	10 g
20% glycerol	50 mL of 100%
0.01% bromophenol blue	0.025 g
125 mM Tris-HCl, pH 6.8	4.91 g

### Washing buffer

<b>Final concentration</b>	<b>For 250 mL</b>
2.5% Triton X-100	6.25 mL of 100%
50 mM Tris-HCl, pH 7.5	12.5 mL of 1 M stock
5 mM CaCl <sub>2</sub>	625 µL of 2 M stock
1 µM ZnCl <sub>2</sub>	2.5 µL of 0.1 M stock

### Incubation buffer

	<b>For 250 mL</b>
Triton x 100 1 %	2.5 mL of 100 %
Tris HCl 50 mM pH 7.5	12.5 mL of 1 M
5mM CaCl <sub>2</sub>	625 µl of 2 M
1 µM ZnCl <sub>2</sub>	2.5 µl of 0.1 M

### Staining solution

	<b>For 100 mL</b>
Methanol	40 mL
Acetic acid	10 mL
H <sub>2</sub> O	50 mL
Coomassie Blue	0.5 g

### Destaining solution

	<b>For 1 L</b>
Methanol	400 mL
Acetic acid	100 mL
H <sub>2</sub> O	500 mL