

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Lysis Buffer Compatibility Table

| LYSIS BUFFER | USAGE | VENDOR AND CATALOG | RANGE TESTED* | CHEM SIGNAL | CHEM RESOLUTION | FLUORESCENT STANDARDS MW SIZING | RECOMMENDATIONS |
|-------------------|---------------------------------|----------------------|---|--|--|---|--|
| T-PER | Tissue Protein | Thermo 78510 | 50–90% | <ul style="list-style-type: none"> 12–230 kDA: Slight signal decrease \geq50% 66–440 kDA: Slight signal decrease \geq50% | <ul style="list-style-type: none"> 2–40 kDA: No effect at 75% 66–440 kDA: Resolution decrease at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: No effect at 75% 66–440 kDA: Signal decrease and smearing at $>$50% | <ul style="list-style-type: none"> 12–230 kDA: For best signal, dilute 1:2 in 0.1X Sample Buffer 66–440 kDA: For best results, dilute 1:3 in 0.1X Sample Buffer |
| M-PER | Mammalian cells | Thermo 78503 | <ul style="list-style-type: none"> 2–40 kDA: 75% 12–230 kDA: 50–900% 66–440 kDA: 50–900% | No effect | No effect | No effect | Samples can be used undiluted with Master Mix |
| RIPA | General Whole Cell Lysis Buffer | Cell Signalling 9806 | 0–90% | <ul style="list-style-type: none"> 2–40 kDA: Increase signal at 75% 12–230 kDA: Signal decrease at 90% 66–440 kDA: Signal loss at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Peak broadening $<$10 kDA at 75% 66–440 kDA: Resolution decreases at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Compression of standards at 75% 12–230 kDA: 230 kDA signal decrease. Non-linear compression causing 1 kDA and 29 kDA standards to run closer 66–440 kDA: Signal decrease and smearing at 50% | <ul style="list-style-type: none"> 2–40 kDA: Prepare Biotinylated Ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards 12–230kDa: For best signal and resolution, dilute 1:2 in 0.1X Sample Buffer 66–440 kDA: Dilute at least 1:3 |
| Cell Lysis Buffer | General Whole Cell Lysis Buffer | Cell Signalling 9803 | 0–90% | <ul style="list-style-type: none"> 2–40 kDA: Increase in signal at 75% 12–230 kDA: Slight signal decrease at high MW region 66–440 kDA: Signal loss at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Peak broadening $<$10 kDA at 75% 66–440 kDA: Resolution decrease at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Compression of standards at 75% 12–230 kDA: 230 kDA signal decrease. Non-linear compression causing 1 kDA and 29 kDA standards to run closer 66–440 kDA: Signal decrease in 280 kDA at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Prepare Biotinylated Ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards 12–230kDa: For best signal and resolution, dilute 1:2 in 0.1X Sample Buffer 66–440 kDA: Dilute at least 1:3 |
| SDS Lysis Buffer | General Whole Cell Lysis Buffer | Millipore #20–163 | 0–90% | No effect | No effect | <ul style="list-style-type: none"> 2–40 kDA: Significant signal decrease for 1 kDA standard and registration due to peak splitting at 75% 12–230 kDA: Smearing of Std 1 peak | <ul style="list-style-type: none"> 12–230 kDA: For best signal and resolution, dilute 1:2 in 0.1X Sample Buffer. Prepare Biotinylated Ladder in same sample buffer for consistent sizing |
| IP Lysis Buffer | General Whole Cell Lysis Buffer | Thermo 87787 | 0–90% | <ul style="list-style-type: none"> 2–40 kDA: Slight signal increase at 75% 12–230 kDA: Slight signal decrease at high MW region 66–440 kDA: Signal decrease at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Peak broadening $<$10 kDA at 75% 66–440 kDA: Resolution decreases at $>$50% | <ul style="list-style-type: none"> 2–40 kDA: Compression of standards at 75% 12–230 kDA: Improved fluorescent standard signal and resolution 66–440 kDA: Signal decrease and smearing in 280 kDA standard at $>$20% | <ul style="list-style-type: none"> 2–40 kDA: Prepare Biotinylated Ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards 12–230 kDA: For best signal and resolution, dilute 1:2 in 0.1X Sample Buffer. Prepare Biotinylated Ladder in same sample buffer for consistent sizing 66–440 kDA: Dilute 1:3 |

*All results compared against ProteinSimple Bicine/CHAPS Buffer.

All results apply to Simple Western size assays with Split Running Buffer (RB) 2 (12–230 kDA), Split RB3 (66–440 kDA), Split RB4 (2–40 kDA) as indicated for each lysis buffer or buffer component tested.

Unless otherwise stated for each separation matrix, the chemi signal, resolution, fluorescent standards, and molecular weight sizing were not affected by the use of addition of the listed components in the ranges tested.

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Lysis Buffer Compatibility Table, continued

| LYSIS BUFFER | USAGE | VENDOR AND CATALOG | RANGE TESTED* | CHEM SIGNAL | CHEM RESOLUTION | FLUORESCENT STANDARDS MW SIZING | RECOMMENDATIONS |
|--------------|---------------------------------|--------------------|---------------|--|--|---|---|
| Cellytic MT | General Whole Cell Lysis Buffer | Sigma C3228 | 0–90% | <ul style="list-style-type: none"> 2–40 kDa: No effect at 75% 66–440 kDa: Signal decrease at 90% | <ul style="list-style-type: none"> 2–40 kDa: No effect at 75% 66–400 kDa: Resolution decrease at 90% | <ul style="list-style-type: none"> 2–40 kDa: No effect at 75% 12–230 kDa: Improved Fluorescent Standard signal and resolution 66–440 kDa: Signal and resolution decrease in 280 kDa at 90% | <ul style="list-style-type: none"> 12–230 kDa: Samples can be used undiluted with Master Mix 66–440 kDa: For best results, dilute 1:2 in 0.1X Sample Buffer |

Lysis Buffer Compatibility Table (for Fluorescence Detection Only)

| LYSIS BUFFER | USAGE | VENDOR AND CATALOG | RANGE TESTED* | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS MW SIZING | RECOMMENDATIONS | PROTEIN NORMALIZATION |
|------------------------|---------------------------------|----------------------|---------------|---|--|--|--|--|
| T-PER | Tissue Protein | Thermo 78510 | 100% | <ul style="list-style-type: none"> 12–230 kDa: Slight signal decrease 66–440 kDa: Slight signal decrease | <ul style="list-style-type: none"> 2–40 kDa: No effect 66–440 kDa: Slight resolution decrease | <ul style="list-style-type: none"> 2–40 kDa: No effect 66–440 kDa: Signal decrease and smearing | | <ul style="list-style-type: none"> 2–40 kDa: No effect 12–230 kDa: No effect 66–440 kDa: No effect |
| RIPA | General Whole Cell Lysis Buffer | Cell Signalling 9806 | 100% | <ul style="list-style-type: none"> 2–40 kDa: Increase in signal at highest concentration of buffer 12–230 kDa: Signal decrease 66–440 kDa: Signal loss | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening such that targets ≤ 10 kDa may be difficult to detect 66–440 kDa: Resolution decrease | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards—1 kDa and 26 kDa run closer 12–230 kDa: 1 kDa and 29 kDa stds run closer 66–440 kDa: Signal decrease and smearing | Prepare biotinylated ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards. | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards causes PN curve to shift but no effect on signal. 12–230 kDa: No effect 66–440 kDa: No effect |
| Cell Lysis Buffer | General Whole Cell Lysis Buffer | Cell Signalling 9803 | 100% | <ul style="list-style-type: none"> 2–40 kDa: Increase in signal at highest concentration of buffer 12–230 kDa: Signal decrease 66–440 kDa: Signal loss | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening such that targets ≤ 10 kDa may be difficult to detect 66–440 kDa: Resolution decrease | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards – 1 kDa and 26 kDa run closer 12–230 kDa: 1 kDa and 29 kDa stds run closer 66–440 kDa: Signal decrease and smearing | Prepare biotinylated ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards. | <ul style="list-style-type: none"> 2–40 kDa: compression of standards causes PN curve to shift but no effect on signal. 12–230 kDa: No effect 66–440 kDa: No effect |
| Pierce IP Lysis Buffer | General Whole Cell Lysis Buffer | Thermo 87787 | 100% | <ul style="list-style-type: none"> 2–40 kDa: Increase in signal at highest concentration of buffer 12–230 kDa: Signal decrease 66–440 kDa: Signal loss | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening such that targets ≤ 10 kDa may be difficult to detect 66–440 kDa: Resolution decrease | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards—1 kDa and 26 kDa run closer 66–440 kDa: Signal decrease and smearing in Std 280 kDa | Prepare biotinylated ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards. | <ul style="list-style-type: none"> 2–40 kDa: compression of standards causes PN curve to shift but no effect on signal. 12–230 kDa: No effect 66–440 kDa: No effect |

*100% of available volume which translates to 76% of absolute volume

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Buffer Reagents

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|--|--|--|--|---|-----------------------|-----------|
| Bicine | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 20–50 mM | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 20–50 mM | No effect | No effect | No effect | No effect |
| TrisCl, pH 7.5 | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 10–50 mM | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 10–50 mM | No effect | No effect | No effect | No effect |
| HEPES, pH 8.0 | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 10–50 mM | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 10–50 mM | 2–40 kDa: Signal increase ≥ 25 mM | 2–40 kDa: Decrease in resolution ≥ 25 mM | No effect | No effect |
| Sodim Phosphate (NaH_2PO_4 / Na_2HPO_4) | <ul style="list-style-type: none"> 2–40 kDa: 7.5–30 mM 12–230 kDa: 10–30 mM | <ul style="list-style-type: none"> 2–40 kDa: 12.5–50 mM 12–230 kDa: 10–50 mM | No effect | No effect | No effect | No effect |
| MES, pH 6.7 | <ul style="list-style-type: none"> 2–40 kDa: 1.25 - 100 mM | <ul style="list-style-type: none"> 2–40 kDa: 1.25 - 100 mM | No effect | No effect | No effect | No effect |

Buffer Reagents (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|---------------|-------------|------------------|---|--|-----------------------|-----------|---|
| HEPES, pH 8.0 | 12.5–50 mM | 12.5–50 mM | <ul style="list-style-type: none"> 2–40 kDa: Increase in signal ≥ 25 mM 12–230 kDa: Signal decrease ≥ 25 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in resolution ≥ 25 mM | No effect | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |
| MES, pH 6.7 | 1.25–100 mM | 1.25–100 mM | No effect | No effect | No effect | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |

Dyes

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|------------------|-------------|------------------|--|--|-----------------------|-----------|
| Bromophenol Blue | 0.001–0.01% | 0.001–0.01% | <ul style="list-style-type: none"> 12–230 kDa: Slight decrease at 0.01% 66–440 kDa: Decrease at low lysate concentration at $>0.1\%$ | <ul style="list-style-type: none"> 66–440 kDa: Decrease in resolution in 0.02 mg/mL HeLa at 0.01% | No effect | No effect |
| Phenol Red | 0.05%–0.01% | 0.05–0.01% | No effect | No effect | No effect | No effect |

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Dyes (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|------------|------------|------------------|--------------------|------------------------|-----------------------|-----------|---|
| Phenol Red | 0.01% | 0.005–0.01% | No effect | No effect | No effect | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |

Salts

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|--------------------|------------|--|---|--|--|---|
| NaCl | 0–700 mM | <ul style="list-style-type: none"> 2–40 kDa: 0–150 mM 12–230 kDa: 0–300 mM 66–440 kDa: 0–300 mM | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening at >150 mM 12–230 kDa: Signal decrease at 300 mM 66–440 kDa: Low signal >300 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in resolution >150 mM 12–230 kDa: Slight decrease at 300 mM 66–440 kDa: Slight decrease at >300 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in signal at >150 mM 12–230 kDa: Std 230 kDa resolution decreases at >300 mM 66–440 kDa: Std 230 kDa slight smear at >700 mM | <ul style="list-style-type: none"> 2–40 kDa: No effect 12–230 kDa: MW sizing may be affected at high MW region due to 230 kDa standard resolution decrease at >300 mM 66–440 kDa: No effect |
| NH ₄ Cl | 0–300 mM | <ul style="list-style-type: none"> 2–40 kDa: 0–100 mM 12–230 kDa: 0–150 mM 66–440 kDa: 0–150 mM | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening at >100 mM 12–230 kDa: Lower signal on high MW end at >150 mM 66–440 kDa: Lower signal on high MW end at >150 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in resolution at >100 mM 12–230 kDa: Loss of resolution >150 mM 66–440 kDa: Loss of resolution >150 mM | <ul style="list-style-type: none"> 12–230 kDa: Std 230 kDa resolution decrease at >75 mM 66–440 kDa: Std 280 kDa resolution decrease at >75 mM | <ul style="list-style-type: none"> 2–40 kDa: No effect 12–230 kDa: MW sizing may be affected at high MW region due to 230 kDa standard resolution decrease at >75 mM 66–440 kDa: No effect |
| MgCl ₂ | 0–10 mM | 0–10 mM | No effect | No effect | No effect | No effect |

Salts (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|--------------------|------------|------------------|---|--|---|-----------|---|
| NaCl | 300 mM | 0–150 mM | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening at >150 mM 12–230 kDa: Signal decrease at 300 mM 66–440 kDa: Low signal >300 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in resolution at >150 mM 12–230 kDa: Slight decrease at 300 mM 66–440 kDa: Slight decrease at 300 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in signal at >150 mM 12–240 kDa: Std 230 kDa resolution decrease 66–440 kDa: Std 280 kDa smearing | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |
| NH ₄ Cl | 150 mM | 0–100 mM | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening at >100 mM 12–230 kDa: Signal decrease at 150 mM 66–440 kDa: Low signal >150 mM | <ul style="list-style-type: none"> 2–40 kDa: Decrease in resolution at >150 mM 12–230 kDa: Resolution decrease at 150 mM 66–440 kDa: Resolution decrease at 300 mM | <ul style="list-style-type: none"> 12–240 kDa: Std 230 kDa resolution decrease 66–440 kDa: Std 280 kDa smearing, resolution decrease | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |
| MgCl ₂ | 10 mM | 0–10 mM | No effect | No effect | No effect | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Reducing Agents

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|-----------|------------|---|--------------|------------------|-----------------------|-----------|
| DTT | 40–80 mM | 40–80 mM | No effect | No effect | No effect | No effect |
| βME | 25–400 mM | 25–400 mM | No effect | No effect | No effect | No effect |
| TCEP | 0.5–2.0mM | <ul style="list-style-type: none"> • 2–40 kDa: 0.5–20 mM • 12–230 kDa: 0.5 mM • 66–440 kDa: 0.5 mM | No effect | No effect | No effect | No effect |

Reducing Agents (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|-----------|------------|------------------|--------------------|------------------------|-----------------------|-----------|-----------------------|
| DTT | 80 mM | 40–80 mM | No effect | No effect | No effect | No effect | No effect |
| BME | 400 mM | 25–400 mM | No effect | No effect | No effect | No effect | No effect |
| TCEP | 20 mM | 0.6–20 mM | No effect | No effect | No effect | No effect | No effect |

Denaturing Agents

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|---------------|--|---|--|--|--|---------------------------|
| Urea | 0–7.6 M | <ul style="list-style-type: none"> • 2–40 kDa: 0–7.6 M • 12–230 kDa: 0–4 M • 66–440 kDa: 0–2 M | 66–440 kDa: Slight decrease at 4 M | 66–440 kDa: Slight decrease at 4 M | 12–230 kDa: Std 1 slight front end smear, but picked up consistently | No effect |
| Urea/Thiourea | <ul style="list-style-type: none"> • 2–40 kDa: 1 M Urea/ 0.2 M Thiourea–5 M Urea/1 M Thiourea • 12–230 kDa: 2 M Urea/0.4 M Thiourea–5 M Urea/1 M Thiourea • 66–440 kDa: 1 M Urea/0.2 M Thiourea–5 M Urea/1 M Thiourea | <ul style="list-style-type: none"> • 2–40 kDa: 1 M Urea/0.2 M Thiourea–5 M Urea/1 M Thiourea • 12–230 kDa: 2 M Urea/0.4 M Thiourea–5 M Urea/1 M Thiourea • 66–440 kDa: 0–2 M Urea/0.4 M Thiourea | 66–440 kDa: Slight decrease at 5 M Urea/1 M Thiourea | 66–440 kDa: Slight decrease at 5 M Urea/1 M Thiourea | 12–230 kDa: Std 1 slight front end smear, but picked up consistently | No effect in range tested |

Denaturing Agents (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|-----------|------------|------------------|--------------------|------------------------|-----------------------|-----------|---------------------------------|
| Urea | 7.6 M | 0–7.6 M | No effect | No effect | No effect | No effect | No effect when FRESHLY prepared |

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Detergents

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|---------------------|--|--|--|--|---|---|
| Triton X-100 | 0–2% | 0–0.5% (All MW ranges) | <ul style="list-style-type: none"> 2–40 kDa: Severe peak broadening at $\geq 1\%$ 12–230 kDa: Signal loss $> 1\%$ 66–440kDa: Signal loss $> 0.5\%$ | <ul style="list-style-type: none"> 2–40 kDa: Slight resolution loss $\geq 1\%$ 12–230 kDa: Slight resolution loss $\geq 1\%$ | <ul style="list-style-type: none"> 2–40 kDa: Peak splitting of 1kDa standard at $\leq 0.25\%$ 12–230 kDa: Std 1 front end smeared, but other standards not affected | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards (1 kDa and 26 kDa run closer) at $> 0.5\%$ 12–230 kDa: Affects sizing at high MW $> 0.5\%$ due to 230 kDa standards resolution loss |
| NP40 | 0–2% | <ul style="list-style-type: none"> 2–40 kDa: 0–0.5% 12–230 kDa: 0–2% 66–440 kDa: 0–2% | 2–40 kDa: Severe peak broadening at $\geq 1\%$ | 2–40 kDa: Poor resolution | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening for Std 1 kDa 12–230 kDa: Slight smear of Std 1, other Stds not affected | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards (1 kDa and 26 kDa run closer) at $> 0.5\%$ 66–440 kDa: Affects sizing $> 2\%$ |
| Igepal CA 630 | <ul style="list-style-type: none"> 2–40 kDa: 0–0.2% 12–230 kDa: 0–1% | <ul style="list-style-type: none"> 2–40 kDa: 0–0.2% 12–230 kDa: 0–1% | No effect | No effect | 2–40 kDa: Peak splitting of 1 kDa standard | 2–40 kDa: May affect MW sizing for small MW targets due to Std 1 splitting |
| C7BZO | <ul style="list-style-type: none"> 2–40 kDa: 0–1% 12–230 kDa: 0–1% 66–440kDa: 0–1% | <ul style="list-style-type: none"> 2–40 kDa: 0–0.025% 12–230 kDa: 0–0.5% 66–440 kDa: 0–0.5% | <ul style="list-style-type: none"> 2–40 kDa: Severe peak broadening at $\geq 0.5\%$ 12–230 kDa: Signal loss $> 1\%$ 66–440 kDa: Signal decreased at 1% for high MW targets | 2–40 kDa: Poor resolution at $\geq 0.5\%$ | 2–40 kDa: Peak broadening for Std 1kDa 12–230 kDa: Slight compression at 1% | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards (1 kDa and 26 kDa run closer) at $> 0.5\%$ 12–230 kDa: May affect sizing at high MW if $> 1\%$ due to compression |
| CHAPS | 0.6–2% | 0.6–2% | No effect | No effect | No effect | <ul style="list-style-type: none"> 2–40 kDa: Compression of standards (1 kDa and 26 kDa run closer) at 2%. Prepare Biotinylated Ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards 12–230 kDa: Prepare Biotinylated Ladder in same sample buffer for accurate MW sizing |
| SDS | 1–2% | 1–2% | 2–40 kDa: Peak broadening at 2% | 2–40 kDa: Decrease resolution at 2% | <ul style="list-style-type: none"> 2–40 kDa: 1kDa Std peak splitting at higher concentrations 12–230 kDa: 1kDa standard slight smear at 1% 66–440 kDa: No effect or slight Std 1 peak splitting) | May affect sizing due to 1 kDa Std peak splitting |
| Sodium Deoxycholate | 0–1% | <ul style="list-style-type: none"> 2–40 kDa: 0–1% 12–230 kDa: 0–0.5% 66–440 kDa: 0–0.5% | 12–230 kDa: Signal decrease $\geq 0.5\%$ | No effect | <ul style="list-style-type: none"> 2–40 kDa: Peak splitting of 1kDa standard at $\geq 0.5\%$ | No effect |
| LDS | <ul style="list-style-type: none"> 2–40 kDa: 0.5–2% 12–230 kDa: 0–1% 66–440 kDa: 0–1% | 0.5–1% | 2–40 kDa: Peak broadening at 2% | 2–40 kDa: Decrease resolution at 2% | <ul style="list-style-type: none"> 2–40 kDa: Peak splitting of 1kDa standard at $\geq 1\%$ 12–230 kDa: 1kDa standard slightly smeared $\geq 1\%$, but consistently picked up | 12–230 kDa: May affect sizing due to Std 1 smear |
| Sarkosyl | 0–1% | <ul style="list-style-type: none"> 2–40 kDa: 0–0.5% 12–230kDa: 0–1% 66–440kDa: 0–1% | 2–40 kDa: Peak broadening at 1% | 2–40 kDa: Decrease resolution at 1% | <ul style="list-style-type: none"> 2–40 kDa: Peak broadening for 1kDa Std peak | No effect in range tested |

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Detergents, continued

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|-----------------|------------|--|---|---|--|--|
| Octyl glucoside | 0–120 mM | <ul style="list-style-type: none"> • 2–40 kDa: Not compatible • 12–230 kDa: 0–60 mM • 66–440 kDa: 0–60 mM | 12–230 kDa: Slight decrease in signal at 120 mM | 12–230 kDa: Slight decrease in resolution at 120 mM | <ul style="list-style-type: none"> • 12–230 kDa: Slight compression at 120 mM | 12–230 kDa: May affect sizing due to compression |

Detergents (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|---------------------|------------|------------------|---|---|---|--|--|
| Triton X–100 | 1% | 0–0.5% | <ul style="list-style-type: none"> • 2–40 kDa: Severe peak broadening at $\geq 1\%$ • 12–230 kDa: Signal loss $\geq 1\%$ • 66–440 kDa: signal loss $\geq 1\%$ | <ul style="list-style-type: none"> • 2–40 kDa: Poor resolution at $\geq 1\%$ • 12–230 kDa: resolution decrease | 2–40 kDa: Peak splitting of 1 kDa standard at $\leq 0.25\%$ | 2–40 kDa: Compression of standards (1 kDa and 26 kDa run closer) at $>0.5\%$ | <ul style="list-style-type: none"> • 2–40 kDa: compression of standards causes PN curve to shift but no effect on signal. • 12–230 kDa: no effect • 66–40 kDa: no effect |
| C7BzO | 1% | 0–0.25% | <ul style="list-style-type: none"> • 2–40 kDa: Severe peak broadening at $\geq 1\%$ • 12–230 kDa: Signal loss $\geq 1\%$ • 66–440 kDa: signal loss $\geq 1\%$ | 2–40 kDa: Poor resolution at $\geq 0.5\%$ | <ul style="list-style-type: none"> • 2–40 kDa: Peak broadening for 1 kDa standard • 12–230 kDa: slight compression at 1% | Compression of standards (1 kDa and 26 kDa run closer) at $\geq 0.5\%$ | <ul style="list-style-type: none"> • 2–40 kDa: compression of standards causes PN curve to shift but no effect on signal. • 12–230 kDa: no effect 6 • 6–40 kDa: no effect |
| CHAPS | 2% | 0.6–2% | No effect | No effect | No effect | 12–230 kDa: Compression of standards (1 kDa and 26 kDa run closer) at 2%. Prepare biotinylated ladder in the same buffer for consistent sizing. Separation time may be increased to offset compression of standards. | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |
| SDS | 2% | 1–2% | <ul style="list-style-type: none"> • 2–40 kDa: Peak broadening at 2% • 66–440 kDa: Signal decrease at 2% | Decreased resolution at 2% | <ul style="list-style-type: none"> • 2–40 kDa: Peak splitting of 1 kDa standard • 12–230 kDa: Std 1 kDa smear slightly | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |
| Sodium deoxycholate | 1% | 0–1% | No effect | No effect | <ul style="list-style-type: none"> • 2–40 kDa: Peak splitting of 1 kDa standard at 1% • 12–230 kDa: Peak splitting of 1 kDa standard at 1% | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |
| LDS | 2% | 0.5–1% | <ul style="list-style-type: none"> • 2–40 kDa: Peak broadening at 2% • 66–440 kDa: Signal decrease at 2% | Poor resolution at 2% | <ul style="list-style-type: none"> • 2–40 kDa: Peak splitting of 1 kDa standard at $\geq 1\%$ • 12–230: std 1 kDa slightly smear • 66–440 kDa: std 230 kDa smearing | No effect | <ul style="list-style-type: none"> • 2–40 kDa: compression of standards causes PN curve to shift but no effect on signal. • 12–230 kDa: no effect • 66–40 kDa: no effect |
| Sarkosyl | 1% | 0–0.5% | 2–40 kDa: Peak broadening at $\geq 1\%$ | 2–40 kDa: Poor resolution at $\geq 1\%$ | <ul style="list-style-type: none"> • 2–40 kDa: Peak broadening for 1 kDa standard • 12–230 kDa: Peak broadening for 1 kDa standard | No effect | No effect in 2–40 kDa, 12–23 kDa and 66–440 kDa |

Simple Western Size Assay Buffer Compatibility (Jess, Abby, Wes, Sally Sue, and Peggy Sue)

Fixative

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|--------------|------------|------------------|--|--|--|-----------|
| Formaldehyde | 0–0.2% | 0–0.1% | 12–230 kDa: Signal decrease $\geq 0.1\%$ | 12–230 kDa: Resolution slightly decrease at $\geq 0.1\%$ | 12–230 kDa: Std 1 front end slightly smeared but picked up consistently $>0.1\%$ | No effect |

Fixative (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|--------------|------------|------------------|--------------------|--|-----------------------|-----------|-----------------------|
| Formaldehyde | 0–0.1% | 0.10% | No effect | 12–230 kDa: Slight resolution decrease | No effect | No effect | No effect |

Viscosity/Density Additives

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|---------------|------------|------------------|---|--|--|--|
| Glycerol | 0–20% | 0–20% | No effect | <ul style="list-style-type: none"> 12–230 kDa: Possible slight decrease in resolution in 0.02 mg/mL HeLa at 20% | 12–230 kDa: Std 1 front end slightly smeared, but picked up consistently | No effect |
| Sucrose | 0–300 mM | 0–300 mM | No effect | No effect | 12–230 kDa: Std 1 front end slightly smeared, but picked up consistently | No effect |
| PEG MW 20,000 | 0–5% | 0–0.05% | <ul style="list-style-type: none"> 2–40 kDa: Extreme peak broadening at $\geq 0.1\%$ 12–230 kDa: Signal loss $\geq 1\%$ 66–440kDa: Signal loss $\geq 1\%$ | Resolution loss $\geq 1\%$ | Complete resolution loss $\geq 1\%$ | Affects sizing due to standards resolution loss. Affect minimized at $\leq 0.05\%$ |

Viscosity/Density Additives (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|---------------|------------|------------------|--|------------------------|---|---------------------------------|-----------------------|
| Glycerol | 20% | 0–20% | No effect | No effect | No effect | No effect | No effect |
| Sucrose | 300 mM | 0–300 mM | No effect | No effect | No effect | No effect | No effect |
| PEG MW 20,000 | 1% | 0–0.06% | <ul style="list-style-type: none"> 2–40 kDa: Extreme peak broadening at $\geq 0.1\%$ | Loss in resolution | <ul style="list-style-type: none"> Extreme peak broadening of 1 kDa standard 12–230 kDa: Peak splitting for 1 kDa std | Minimal effect at $\leq 0.06\%$ | No effect |

Miscellaneous Reagents

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|-----------|------------|---|--|------------------|--|-----------|
| EDTA | 0–40 mM | 0–40 mM | No effect | No effect | No effect | No effect |
| Imidazole | 0–100 mM | <ul style="list-style-type: none"> • 2–40 kDa: 0–100 mM • 12–230 kDa: 0–50 mM • 66–440 kDa: 0–100 mM | <ul style="list-style-type: none"> • 12–230 kDa: Signal decrease at 100 mM especially for high MW targets | No effect | <ul style="list-style-type: none"> • 12–230 kDa: Std 1 front end slightly smeared, but picked up consistently | No effect |

Miscellaneous Reagents (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|-----------|------------|------------------|----------------------------------|------------------------|-----------------------|-----------|-----------------------|
| EDTA | 40 mM | 0–40 mM | 2–40 kDa: Slight signal increase | No effect | No effect | No effect | No effect |

Buffer

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | CHEMI SIGNAL | CHEMI RESOLUTION | FLUORESCENT STANDARDS | MW SIZING |
|-----------|------------|------------------|---|--|--|--|
| PBS | 50–90% | 50% | <ul style="list-style-type: none"> • 12–230 kDa: Slight loss of signal at 90% • 66–440kDa: Slight loss of signal at 90% | <ul style="list-style-type: none"> • 12–230 kDa: Slight loss of resolution at 90% | <ul style="list-style-type: none"> • 12–230 kDa: Slight smearing of Std 230 kDa at 90% • 66–440 kDa: Slight smearing of Std 280 kDa at 90% | <ul style="list-style-type: none"> • 12–230 kDa: May affect sizing at HMW at 90%, due to smearing of Standard 230 kDa |

Buffer (for Fluorescence Detection Only)

| COMPONENT | TEST RANGE | COMPATIBLE RANGE | FLUORESCENT SIGNAL | FLUORESCENT RESOLUTION | FLUORESCENT STANDARDS | MW SIZING | PROTEIN NORMALIZATION |
|-----------|------------|------------------|------------------------------------|---|--|-----------|-----------------------|
| PBS | 1X | 0.25X–1X | 66–440 kDa: Slight signal decrease | <ul style="list-style-type: none"> • 12–230 kDa: Slight loss of resolution | <ul style="list-style-type: none"> • 66–440kDa: Std 280 kDa smearing slightly | No effect | No effect |



Toll-free: (888) 607-9692
Tel: (408) 510-5500
info@proteinsimple.com
proteinsimple.com

© 2021 ProteinSimple. ProteinSimple, Simple Western, Jess, Abby, Wes, Peggy Sue, Sally Sue and the ProteinSimple logo are trademarks and/or registered trademarks of ProteinSimple.