

# 10 Tips for Top Checkweigher Performance

Setting up your checkweigher to achieve optimum performance isn't a luxury—it is critical to meeting your production obligations. This means knowing the signs of when and where you need to take action—and what action to take quickly and effectively to get back in the zone of peak operation. Minimize give-away while making sure all your products meet labeled weight and nutritional information. Don't be a bottleneck!



## 1. Mechanical Alignment

Is the frame firmly on the ground? Are the conveyors all level and aligned? Without correct mechanical alignment the weighing process will fail to meet specifications.



## 2. Photoeyes and Rejectors

All photoeyes should be checked for position and operation. Rejectors should fire at the proper time to ensure the rejected product goes into the reject bin properly.



## 3. Belts or Chains

Inspect your belts or chains for wear or product build-up. Replace if needed. Otherwise inaccurate weights are guaranteed. Inspect timing belts for wear, cracks or missing teeth.



## 4. Product Transfer

Do products jump or vibrate when they transfer onto the checkweigher? Are the gaps between the conveyors minimized? Stable products weigh better! Are the infeed and weigh table belts running at the same speeds? This is critical.



## 5. Product Pitch

Is there enough gap between products to guarantee there is never more than one product on the weigh table at a time during the weighing process? You may need to speed up the infeed to correct the issue.



## 6. Product Setup

Make sure the proper dimensions and weight of the product are correctly entered into the product recipe. Are the cutpoints for over-weight and under-weight products set to correct product specifications?



## 7. Calibration and Auto Zero

Has the checkweigher been calibrated recently for zero and span both empty and with the product? A calibration should be done daily or whenever switching product. Is there enough gap during production for the auto zero routine to adjust?



## 8. Grey Zone Validation

When was your calibration last checked? A grey zone check should be run before every shift or product change-over to ensure optimum weights every day. Be sure to run a package at least 20 times to get a statistically valid result.



## 9. Alarms

Are the triggers for alarm conditions set properly? An alarm that constantly goes on will be ignored by operators.



## 10. Reports

Reports that deliver lot, batch or shift information on accept/reject quantities need to be formatted properly to be accurate representations of production.