

# Aerospace aluminum alloy 2099 friction stir weld analysis

GKN Aerospace utilizes transmission electron microscopy to understand the processing-structure-property relationship

## Introduction

Friction stir welding (FSW) is a joining technique that creates a weld by stirring solid metal between adjoining plates rather than melting it. This solid-state process will locally heat the metal to around the hot forging temperature, which is much lower than traditional fusion welding processes. FSW is broadly utilized in aluminum alloys as it achieves better mechanical properties in comparison to fusion based processes. Yet welding precipitation-hardened aluminum alloys remains a challenge. Using FSW for such alloys results in a reduction on the potency of solidification cracks, lower residual stresses and part distortion. Typically, the process is done via using conventional (CNC) machine tools which is preferential in terms of the process flow and industrialization.

The FSW demonstrations were made at TWI and EWI following guidelines from GKN Aerospace. Replacing drilling and riveting by continuous FSW removes all the stress concentration points due to the occurrence of the drilled holes. It also removes the potential for corrosion due to the presence of a different material (rivet) with respect to the aluminum part.

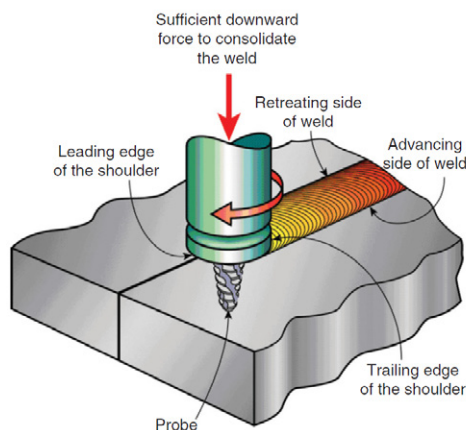


Figure 1: Single-sided stir friction welding site. Image courtesy of TWI.

Multiple weld tests were conducted with variables such as single- or double-sided FSW, machine speed and metal surface treatment. Test samples were provided to Warwick Manufacturing Group (WMG) for analysis by optical microscopy, electron backscattered diffraction (EBSD), and transmission electron microscopy, though a complete understanding remained elusive.

## Methods and results

Hardness mapping revealed to GKN Aerospace the impact of FSW on mechanical performance, where the weld region suffered about 18% loss of harness and strength.

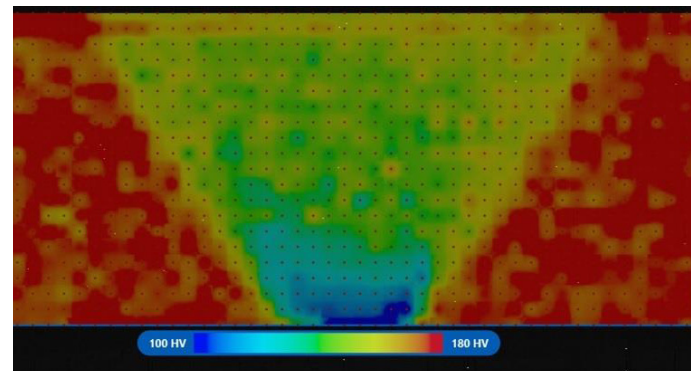


Figure 2. Hardness test mapping on a FSW cross section. Courtesy GKN Aerospace.

WMG performed EBSD imaging using the Versa 3D DualBeam, formerly produced by FEI, which helped provide a deeper dive into the microscopic structure showing detailed differences in grain size in the different regions. GKN Aerospace could then see that the grain structure of the weld had been entirely changed.

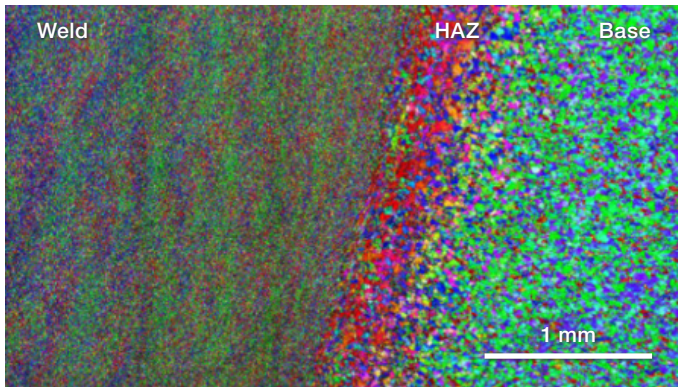


Figure 3. EBSD imaging of the weld, HAZ, and base metal regions, courtesy of WMG (Warwick Manufacturing Group), University of Warwick.

The stirred zone revealed a banded texture with grains finer than the base metal. The HAZ yielded a new grain orientation with some grain growth, and the base metal remained unaffected by the welding process.

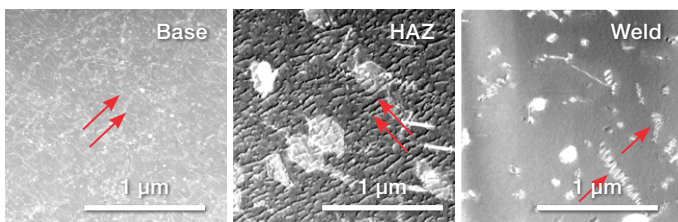


Figure 4: Thermo Scientific Talos F200X STEM mode imaging of dislocations.

The high misorientations of the base metal were indications of plastic deformation, work-hardening, and accumulation of dislocations. These characteristics, which were responsible for the high strength of the Al 2099 base metal, were now absent within the stirred zone. The apparent loss of dislocation substructures and precipitates in the stirred zone negatively affects the strength and ductility.

The Thermo Scientific™ Helios™ 5 Laser PFIB was used to prepare TEM lamellae for nanometer-scale analysis.

TEM analysis using the Thermo Scientific Talos™ F200X TEM imaging in STEM mode across the different regions showed the change in dislocations from base to HAZ to weld.

GKN Aerospace concluded that the base metal and HAZ have a high density of edge dislocations, whereas the weld has lost these in favor of only a few screw dislocations.

Finally, chemical mapping by EDS and particle quantification by Automated Particle Workflow (APW) was conducted. Both the base metal and HAZ showed the  $T_1$  ( $Al_2CuLi$ ) type of precipitate, though they had become coarser in the HAZ. Conversely, the weld area had none of the  $T_1$  precipitates, as the copper is now present in the  $T_\beta$  ( $Al_{7.5}Cu_4Li$ ) precipitate.

If the base metal was considered the optimum structure, then the HAZ retained some of the strengthening  $T_1$  precipitates, and the weld had lost the  $T_1$  precipitate altogether.

### Conclusions

Friction stir welding of aluminum alloys remains a challenging process and there is significant potential for improvement. Optimizing the post-weld heat treatments could provide some of the strength which is lost in the weldment. High-resolution TEM/EDS analysis of precipitates was made simple with the Talos F200X TEM. The Automated Particle Workflow (APW) characterized the change in precipitate size and morphology across the base metal, heat affected zone and weld.

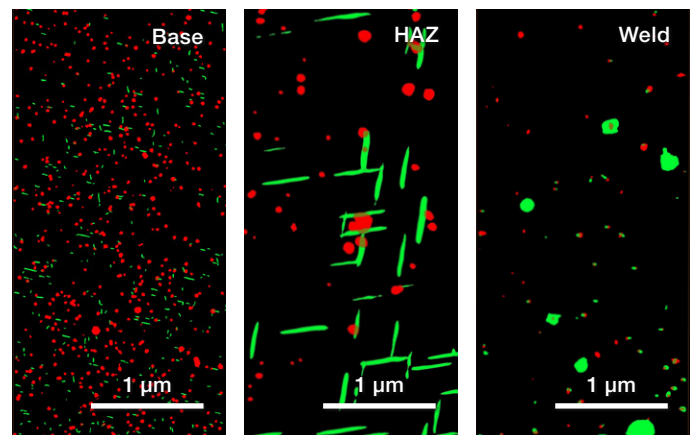



Figure 5: Friction stir welded Al 2099. The green represents copper, the red represents zirconium, and the green platelets are the  $T_1$  precipitate in the base and HAZ.



## Automated Particle Workflow on Aluminum 2099 alloy

Duration 0:53

Find out more at [thermofisher.com/metals-research](https://thermofisher.com/metals-research)

**ThermoFisher**  
SCIENTIFIC

For research use only. Not for use in diagnostic procedures. For current certifications, visit [thermofisher.com/certifications](https://thermofisher.com/certifications)

© 2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AN0176-EN-12-2021